



User Manual

PrecisION[®] Ion Channel Cell Lines

Validated Cell Lines for Use in Cell-Based Functional Assays Including Manual and Automated Patch Clamp or Fluorescent Platforms



Document Number 70-430 Revision 2.0



Table of Contents

Important: Please read this entire user manual before proceeding with the assay.

Overview	1
PrecisION Cell Lines.....	1
Materials Provided.....	1
Storage Conditions	1
Additional Materials Required	1
Unpacking Cell Cryovials.....	2
Cell Culture Protocol.....	3
Cell Thawing	3
Cell Recovery	4
Cell Propagation	5
Cell Cryopreservation	5
Recommendations for Testing	7
1. Preparation of Cells	7
2. Compound Preparation	7
Appendix 1: Complete Media Formulations – Parental Cell Lines.....	8
Appendix 1: Complete Media Formulations – Recombinant Cell Lines	9
Appendix 2: Cell Line Specific Seeding Guidelines.....	15
Document Revision History.....	18
Limited Use License Agreement	19

For additional information or Technical Support, see contact information at the bottom of this page.

Overview

PrecisION cell lines are designed to provide pharmacologically relevant cell models for research and screening applications.

PrecisION Cell Lines

PrecisION cell lines stably express the target of interest in cell backgrounds that are highly suitable for electrophysiology applications.

Materials Provided

Components	
2 vials of cells	Refer to cell line-specific datasheet for shipped cell density

Storage Conditions

Cells are shipped on dry ice and should arrive in a frozen state. To ensure maximum cell viability, store the vials of cells in the vapor phase of liquid nitrogen as soon as possible upon receipt. Please contact technical support immediately, if the cells received were already thawed.



Contact technical support immediately if cells received were already thawed.

- Short-term (less than 24 hours): Store vials at -80°C immediately upon arrival. (DO NOT store at -80°C for longer than 24 hours).
- Long-term (longer than 24 hours): Vials should ONLY be stored in liquid nitrogen (vapor phase ONLY) or in ultra-low temperature (-150°C) freezers.

Additional Materials Required

Refer to the table below to determine the appropriate media and reagents required for the specific cell line used in the assay. The cell type is provided in the cell line-specific datasheet.

Material	Ordering Information Recommended Products (or Equivalent)
Cell line-specific Complete Media	Refer to the tables in Appendix 1 for cell line-specific media requirements
PBS, pH 7.4 (without calcium and magnesium)	ThermoFisher 10010-031 (or equivalent)
AssayComplete™ Cell Detachment Reagent	Eurofins DiscoverX Cat # 92-0009
Single and multichannel micropipettes and pipette tips	
50 mL and 15 mL polypropylene tubes	
1.5 mL microtubes	
Tissue culture disposable pipettes (1 mL-25 mL) and tissue culture flasks (T25, T75 and T225 flasks, etc.)	
Cryovials for freezing cells	
Automated Cell Counter (or Hemocytometer)	
Humidified tissue culture incubator (37°C and 5% CO ₂)	

Unpacking Cell Cryovials

Two cryovials are shipped on dry ice, and each contain cells in 1 mL of Freezing Reagent (refer to the cell line-specific datasheet for the number of cells provided in the vials). The following procedures are for safe storage, handling and removal of cryovials from the vapor phase of liquid nitrogen storage.

1. Cells must arrive in a frozen state on dry ice.



Contact technical support immediately if cells received were already thawed.

2. Frozen cells must be transferred to either the vapor phase of liquid nitrogen or a -80°C freezer immediately upon arrival. If cells will be thawed and used within 24 hours, they can be stored temporarily at -80°C. For long-term storage, cryovials must be stored in the vapor-phase of liquid nitrogen.



Cryovials should be stored in the vapor phase of liquid nitrogen. They are not rated for storage in the liquid phase.

3. Using tongs, remove the cryovials from the liquid nitrogen storage, and place them immediately in dry ice, in a covered container.



A face shield, gloves, and a lab coat should be worn during these procedures.

4. Wait for at least one minute for any liquid nitrogen that may be present inside the vial to evaporate.
5. Proceed with the thawing and propagation protocol in the following section. Refer to the cell line-specific datasheet for appropriate products mentioned in the following protocols.

Cell Culture Protocol

The following procedures are for thawing adherent cells from cryovials, seeding and expanding the cells, and freezing the cells once they have been propagated. Refer to the cell-line specific datasheet to identify the cell type of the cell line. Refer to [Appendix 1](#) to determine the correct complete media formulation for the specific cell type.

Cell Thawing

The following is a protocol for thawing cells into a T75* flask. (*T25 flask where noted)

1. Pre-warm the complete media in a 37°C water bath for 15 minutes, as indicated on the cell line-specific datasheet.
2. Add 9 mL of the complete media into a 15 mL conical tube inside a sterile tissue culture hood. Set aside for [Step 6](#).
3. Remove the cryovials from -80°C or liquid nitrogen vapor storage and immediately place them in dry ice prior to thawing.



Safety Warning: A face shield, gloves, and a lab coat should be worn at all times when handling frozen vials. When removing cryovials from the vapor phase of liquid nitrogen storage, use tongs to place them immediately in dry ice in a covered container. Wait at least 1 minute for any liquid nitrogen inside the vial to evaporate. Do not touch the bottom of the vials at any time, to avoid inadvertent thawing of the cells.

4. Place the cryovials containing the cells in a 37°C water bath briefly, until only small ice crystals remain and the cell pellet is almost completely thawed. The thawing time typically ranges from 1.5 to 2.5 minutes. DO NOT vortex freshly thawed cells.



Do not leave the frozen cell vials unattended in the water bath. Prolonged thawing at 37°C may result in cell death.

5. Decontaminate the external surface of the vials by spraying and wiping them with 70% ethanol. Transfer the vials to a tissue culture hood.
6. Using a sterile pipette, gently transfer the thawed cells to the pre-filled 15 mL conical tube.
7. Rinse the vial with 1 mL of complete media to maximize cell recovery and pipette it into the 15 mL conical tube.
8. Centrifuge the 15 mL conical tube containing cells at 250 x g for 4 minutes.
9. Add 18 mL of complete media to a T75 flask (or 6 mL to a T25 flask for the cell lines noted below). Some cell lines recover better when thawed into a T25 flask (CYL3054, CYL3056, and CYL3096).
10. After centrifugation, aspirate the media from the conical tube with care not to disturb the cell pellet.
11. Gently resuspend the pellet with 2 mL of complete media.
12. Transfer the cell suspension to the T75 (or T25 where noted) flask and gently swirl the flask to distribute the cells evenly in solution.
13. Incubate the flask in a 37°C and 5% CO₂ humidified cell culture incubator.
14. Maintain the cells in culture until they are 60-75% confluent in the T75 (or T25 flask). Do not exceed 80% confluency, as target down-regulation may occur.
15. When passaging cells, refer to the instructions in the [Cell Recovery](#) section.

16. We recommend maintaining the cultures in T225 flasks, being careful not to exceed 80% confluency. All seeding densities in this manual are those recommended for T225 flasks. If other sized flasks are used, adjust seeding densities appropriately, according to the surface area of the flask.

Cell Recovery

The following is a protocol for ensuring maximal cell recovery once the cells become 60-80% confluent in a T225 flask.

1. Pre-warm the complete media in a 37°C water bath for 15 minutes.
2. Remove the T225 flask from the tissue culture incubator and place it in a sterile tissue culture hood.
3. Gently aspirate the media from the T225 flask.
4. Add 15 mL of PBS into the T225 flask, and gently rock the flask back and forth to rinse the cells.
5. Gently aspirate PBS from the flask.
6. Add 4 mL of pre-warmed Cell Detachment Reagent to the flask. Gently rock the flask back and forth to ensure that the inner surface of the flask is uniformly covered with Cell Detachment Reagent.
7. Incubate the flask at 37°C and 5% CO₂ until the cells have detached. If the cell line is a HEK 293 background, the maximum amount of time for Cell Detachment Reagent to be on the cells is about 2 minutes. If the cell line is a CHO-K1 background, the maximum amount of time is about 3 minutes.



The maximum amount of time for Cell Detachment Reagent to be on the cells is different depending on the background. Do not leave Cell Detachment Reagent on cells longer than necessary.

8. Remove the flask from the incubator and observe the cells under a microscope to confirm cell detachment. If necessary, gently tap the edge of the flask to detach cells from the surface. If most of the cells have not yet detached, then return the flask to the incubator for additional 1 minute. Repeat this step until all cells are in suspension.



Some cell lines are highly adherent and therefore may require additional dissociation time.

9. Add 10 mL of pre-warmed complete media to a 50 ml centrifuge tube.
10. Using a pipette, slowly add 16 mL of the complete media to the detached cells in T225 flask to gently rinse them into the suspension. Gently pipet the cell suspension up and down several times to generate a single cell suspension without any clumps.
11. Count the number of viable cells using an automated cell counter or hemocytometer. Determine the number of cells and cell suspension volume required to seed into a T225 flask. Cells should be initially seeded according to [Appendix 2](#). Please Note: These are initial seeding recommendations for the first passages out of thaw. Seeding densities should be adjusted based on the desired time window for expression, generally recommended to be 48 - 72 hours. Cell growth rates will vary depending on the amount of time in continuous culture, and the density of cells prior to lifting them during recent passages.
12. Add 15 mL of growth media to a new T225 flask, followed by addition of the cell suspension. Add an additional volume of the growth media to reach a final volume of 54 mL for a T225 flask.
13. Transfer the flask to a tissue culture incubator and incubate the cells at 37°C and 5% CO₂.

Cell Propagation

The cells can be propagated after a successful recovery, which can be determined by monitoring them under a microscope. Healthy cells should adhere uniformly to the surface of the flask, with only a few cells remaining in suspension.




To maintain logarithmic growth of the cells, cultures should be maintained in a sub-confluent monolayer.

1. If the cells do not appear to be healthy, or if confluency is <25% (Ion Channel cell lines may demonstrate unusual morphology at low densities and within 72 hours of thawing):
 - 1.1. Incubate the flask for additional 24 to 48 hours to allow for cell recovery before exchanging the cell culture media. If culture has been static for more than 3 days, a media exchange may be required. Some lines may require up to 5 days to recover from being thawed.
2. Once the cells have reached 65-75% confluency, split the cells every 2 to 3 days, based on the doubling time of the cell line. Refer to [Appendix 2](#) for cell seeding recommendations.

Cell Cryopreservation

The following procedure is for freezing cells that have been propagated in T225 flasks. This protocol assumes that the cells have reached 65-75% confluency in a sufficient number of flasks to fill the required number of cryovials. The cells will be resuspended in ice-cold CryoMedium (described in [Step 15](#)), and will be cryopreserved in aliquots of 1 mL/vial, at the desired number of cells per vial (e.g. 2.0×10^6 per vial). Refer to the table below for the ratio of components used to prepare the CryoMedium:

Basal Media (Without Selection Antibiotics)	40%
Fetal Bovine Serum (FBS)	50%
Cell-culture grade Dimethyl Sulfoxide (DMSO)	10%

1. Remove T225 flasks from the incubator and place them in a sterile tissue culture hood.
 -  Care should be taken while handling flasks to avoid contamination.
 2. Slowly aspirate the media from the flasks.
3. Add 15 mL of PBS into each T225 flask, and gently rock the flask back and forth to ensure that the cells are rinsed.
4. Gently aspirate PBS from the flask.
5. Add 4 mL of pre-warmed Cell Detachment Reagent to T225 flask.
6. Gently rock the flask back and forth to ensure that the inner surface of the flask is uniformly covered with Cell Detachment Reagent.
7. Incubate the flask at 37°C and 5% CO₂ until the cells have detached. If the cell line is a HEK 293 background, the maximum amount of time for Cell Detachment Reagent to be on the cells is 2 minutes. If the cell line is a CHO-K1 background, the maximum amount of time is 3 minutes.



The maximum amount of time for Cell Detachment Reagent to be on the cells is different depending on the background. Do not leave Cell Detachment Reagent on cells longer than necessary. Doing so may disrupt the ion channel function.

8. Remove the flask from the incubator and view under a microscope to confirm that the cells have detached. If necessary, gently tap the edge of the flask to detach cells from the surface. If most of the cells have not

detached, then return the flask to the incubator for additional 1 minute and repeat this step until all cells are in suspension.

9. Add 16 mL of the cell culture media to each T225 flask to neutralize the Cell Detachment Reagent
10. Gently pipet the cell suspension up and down several times to generate a single cell suspension without any clumps.
11. Transfer the cell suspension from the T225 flask into a 50 mL conical centrifuge tube. If necessary, add an additional volume of the cell culture media to the flasks and rinse to collect the remaining cells, then transfer it into the conical tube. Slowly pipette up and down several times to ensure that a single cell suspension is formed.
12. To determine the number of cells in the suspension:
 - 12.1. Set aside a small fraction (0.5 mL or less) of the suspended cells in a separate tube.
 - 12.2. Transfer an appropriate portion of this fraction to a cell counter or hemocytometer.
 - 12.3. Count cells, calculate the concentration of cells in the suspension, and then calculate the total number of cells in the 50 mL centrifuge tube.
13. Centrifuge the collected cells at 200 x g for 4 minutes.
14. After centrifugation, discard the supernatant, being careful not to disturb the cell pellet.
15. Based on the total cell number calculated in **Step 12**, resuspend the cells to the desired concentration (e.g. 2.0×10^6 cells/mL) with ice-cold CryoMedium.



Keep cells on ice during this process to protect cell viability.

16. Aliquot 1 mL of the cell suspension into each of the labeled 2 mL cryovials. Seal the cryovials tightly.
17. Freeze cells in a -80°C freezer at a controlled rate of -1°C/minute overnight in a dedicated cell freezer or commercially-available freezing chamber. For short-term storage, vials can be stored in the -80°C freezer for a maximum of two weeks.
18. Transfer the vials into the vapor phase of liquid nitrogen for long-term storage.

Recommendations for Testing

General guidelines are provided for cell seeding but may vary based on recording conditions and equipment used for assay. For running the assay, refer to relevant protocols as recommended by the equipment manufacturer or published literature.

1. Preparation of Cells

The following protocol provides steps for harvesting and preparing adherent cells. This protocol assumes that cells have reached a 65-75% confluency and have been cultured in the recommended cell culture media. For cell line-specific information on the control compound, incubation times and temperature schedule, refer to the cell line-specific datasheet.

- 1.1. Ensure that the cells are in the logarithmic growth phase prior to seeding for use in the assay.
- 1.2. Warm the cell culture media in a clean 37°C water bath for 15 minutes.
- 1.3. Gently harvest the cells with the method-specific dissociation and assay preparation protocol.

2. Compound Preparation

For compounds prepared in DMSO, ensure final DMSO concentration in assay is < 0.3 – 0.5%.

Appendix 1: Complete Media Formulations – Parental Cell Lines

Growth media used for the parental cell lines are listed in the following table.

Complete Media	Target	Catalog Number	Component	Supplier Name	Supplier Part Number
HEK Wild Type Media	HEK 293 WT1	CYL3001	DMEM/F-12 (with L--Glutamine)	ThermoFisher	10565-018
	HEK 293 WT2	CYL3002	10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
CHO Wild Type Media	CHO-K1 WT	CYL3000	F-12 (Ham's) + L--Glutamine	ThermoFisher	3176-035
			10% FBS	Millipore Sigma	F2442

Appendix 1: Complete Media Formulations – Recombinant Cell Lines

Growth media used for each cell line background are listed in the following tables.

Complete Media	Target	Catalog Number	Component	Supplier Name	Supplier Part Number
HEK Media 1	hNav1.5	CYL3004	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
	hNav1.1	CYL3009			
	hNav1.6	CYL3010	10% FBS	Millipore Sigma	F2442
	hNav1.7	CYL3011	1% Non-Essential Amino Acids	ThermoFisher	11140-050
	hNav1.4	CYL3024			
	hKir2.1	CYL3032	400 µg/mL G418	Eurofins DiscoverX	92-0030
	hERG-HEK	CYL3039			
	hHCN1	CYL3040			
	hHCN2	CYL3041			
	hHCN3	CYL3042			
	hGluR6	CYL3049			
	hASIC3	CYL3055			
	hTRPV1	CYL3063			
	hTRPV4	CYL3064			
	hTRPV3	CYL3065			
	hTRPA1	CYL3066			
	hCav3.2	CYL3075			
	hCFTR	CYL3088			
	hKv12.2	CYL3089			

HEK Media 2	hNav1.8/ β1	CYL3025	DMEM/F-12 (with L--Glutamine)	ThermoFisher	10565-018
	hGABAA α1/β3/γ2	CYL3053	10% FBS	Millipore Sigma	F2442
		CYL3054 (Thaw into T25 flask)	1% Non-Essential Amino Acids	ThermoFisher	11140-050
	CYL3068		400 µg/mL G418	Eurofins DiscoverX	92-0030
	hGABAA α3/β3/γ2	CYL3073	100 µg/mL Hygromycin	Eurofins DiscoverX	92-0029
	hGABAA α5/β3/γ2		CYL3085	0.625 µg/mL Puromycin	Eurofins DiscoverX
	hGABAA α4/β3/γ2	CYL3086			
	hnAChR α4/α6/β2	CYL3107			
HEK Media 3	hKv7.4	CYL3092	DMEM/F-12 (w/ L--Glutamine)	ThermoFisher	10565-018
	hKv7.4/K v7.5	CYL3096 (Thaw into T25 flask)	10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
	hnAChR α4/β2	CYL3106	200 µg/mL Hygromycin	Eurofins DiscoverX	92-0029
HEK Media 4	hnAChR α7/ric3	CYL3097	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
			10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 µg/mL G418	Eurofins DiscoverX	92-0030
			0.625 µg/mL Puromycin	Eurofins DiscoverX	92-0028

HEK Media 5	nAChR $\alpha 1/\beta 1/\delta/\epsilon$	CYL3052	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
	hnAChR $\alpha 3/\beta 4$	CYL3057	10% FBS	Millipore Sigma	F2442
	hKv4.3/K ChiP2	CYL3069	1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 μ L/mL G418	Eurofins DiscoverX	92-0030
			100 μ g/mL Hygromycin	Eurofins DiscoverX	92-0029
HEK Media 6	hKv7.3/K v7.5	CYL3060	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
	hKir6.2/S UR2A (KATP)	CYL3099	10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 μ g/mL G418	Eurofins DiscoverX	92-0030
			200 μ g/mL Hygromycin	Eurofins DiscoverX	92-0029
HEK Media 7	hGlyRA1	CYL3056 (Thaw into T25 flask)	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
			10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 μ g/mL G418	Eurofins DiscoverX	92-0030
			30 μ M Picrotin	Millipore Sigma	Y0001508

HEK Media 8	Cav1.2 α1C/β2a/ α2δ1	CYL3051	DMEM/F-12 (with L-Glutamine)	ThermoFisher	10565-018
			10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 µg/mL G418	Eurofins DiscoverX	92-0030
			100 µg/mL Hygromycin	Eurofins DiscoverX	92-0029
			0.625 µg/mL Puromycin	Eurofins DiscoverX	92-0028
			144 nM Nitrendipine	Millipore Sigma	N144
CHO Media 1	hNav1.3 hHCN4 hNav1.2	CYL3003 CYL3012 CYL3023	IMDM (with L-Glutamine)	ThermoFisher	12440-053
			10% Dialyzed FBS	ThermoFisher	26400-044
			1% HT Supplement	ThermoFisher	11067-030
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 µg/mL G418	Eurofins DiscoverX	92-0030

CHO Media 2	hKv1.1	CYL3014	F-12 (Ham's) + L-Glutamine	ThermoFisher	31765-035
	hKv1.2	CYL3015	10% FBS	Millipore Sigma	F2442
	hKv1.3	CYL3016			
	hKv1.4	CYL3017	400 µg/mL G418	Eurofins DiscoverX	92-0030
	hKv1.5	CYL3018			
	hKv1.6	CYL3019			
	hKv1.7	CYL3020			
	hKv1.8	CYL3021			
	hKv2.1	CYL3022			
	hERG- CHO	CYL3038			
	hKv3.1	CYL3043			
	hKv3.2	CYL3044			
	hKv3.3	CYL3045			
	CHO Media 3	hKCNQ1 /minK	CYL3007	IMDM (with L-Glutamine)	ThermoFisher
10% Dialyzed FBS				ThermoFisher	26400-044
1% HT Supplement				ThermoFisher	11067-030
1% Non-Essential Amino Acids				ThermoFisher	11140-050
400 µg/mL G418				Eurofins DiscoverX	92-0030
100 µg/mL Hygromycin				Eurofins DiscoverX	92-0029

CHO Media 4	hKv4.2/K CHiP2	CYL3026	F-12 (Ham's) + L-Glutamine	ThermoFisher	31765-035
	hKv7.2/K v7.3	CYL3059	10% FBS	Millipore Sigma	F2442
	hKv2.1/K v9.2	CYL3067	400 µg/mL G418	Eurofins DiscoverX	92-0030
			100 µg/mL Hygromycin (Gibco)	Eurofins DiscoverX	92-0029
CHO Media 5	hKv4.3/K CHiP1	CYL3027	DMEM (with L-Glutamine)	ThermoFisher	11966-025
			10% FBS	Millipore Sigma	F2442
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 µg/mL G418	Eurofins DiscoverX	92-0030
ND7-23 Media	rNav1.8	CYL3050	DMEM (with L-Glutamine)	ThermoFisher	11966-025
			10% Heat-Inactivated FBS**	Millipore Sigma	F2442
			**Heat inactivation must be performed by user by incubating in a 60° water bath for at least 30 minutes.		
			1% Non-Essential Amino Acids	ThermoFisher	11140-050
			400 µg/mL G418	Eurofins DiscoverX	92-0030

Appendix 2: Cell Line Specific Seeding Guidelines

Suggested seeding guidelines for each cell line background are listed in the following tables. Growth rate may vary due to amount of time in culture, and percent confluency at recent passages.

Catalog Number	Cell Line	Seeding Density (x 10⁶) T225 flasks (48-72 hours)
CYL3000	CHO-K1 WT	6.0
CYL3001	HEK WT1	1.8
CYL3002	HEK WT2	1.8
CYL3003	hNav1.3	2.6
CYL3004	hNav1.5	5.3
CYL3007	hKCNQ1/minK	1.8
CYL3009	hNav1.1	5.3
CYL3010	hNav1.6	5.3
CYL3011	hNav1.7	5.3
CYL3012	hHCN4	2.0
CYL3014	hKv1.1	2.0
CYL3015	hKv1.2	2.0
CYL3016	hKv1.3	2.1
CYL3017	hKv1.4	2.1
CYL3018	hKv1.5	2.1
CYL3019	hKv1.6	2.1
CYL3020	hKv1.7	2.1
CYL3021	hKv1.8	2.1
CYL3022	hKv2.1	4.2

CYL3023	hNav1.2	2.6
CYL3024	hNav1.4	4.2
CYL3025	hNav1.8/β1	2.4
CYL3026	hKv4.2/KChIP2	2.1
CYL3027	hKv4.3/KChIP1	2.1
CYL3032	hKir2.1	2.1
CYL3038	hERG	2.1
CYL3039	hERG	5.3
CYL3040	hHCN1	8.3
CYL3041	hHCN2	8.3
CYL3042	hHCN3	2.0
CYL3043	hKv3.1	2.1
CYL3044	hKv3.2	2.1
CYL3045	hKv3.3	2.1
CYL3049	hGluR6	6.8
CYL3050	rNav1.8	2.4
CYL3051	Cav1.2 α1C/β2a/α2δ1	4.2
CYL3052	nAChR α1/β1/δ/ε	2.4
CYL3053	hGABAA α1/β3/γ2	6.8
CYL3054	hCav2.2 α1B/β3/α2δ1	6.0
CYL3055	hASIC3	2.4
CYL3056	hGlyRA1	6.8

CYL3057	hnAChR α 3/ β 4	6.0
CYL3059	hKv7.2/Kv7.3	2.1
CYL3060	hKv7.3/Kv7.5	2.1
CYL3063	hTRPV1	5.4
CYL3064	hTRPV4	2.4
CYL3065	hTRPV3	2.4
CYL3066	hTRPA1	2.4
CYL3067	hKv2.1/Kv9.2	2.1
CYL3068	hGABAA α 3/ β 3/ γ 2	2.3
CYL3069	hKv4.3/KChiP2	2.1
CYL3073	hGABAA α 5/ β 3/ γ 2	2.3
CYL3075	hCav3.2	1.8
CYL3085	hGABAA α 4/ β 3/ γ 2	3.0
CYL3086	hGABAA α 6/ β 3/ γ 2	4.5
CYL3088	hCFTR	1.8
CYL3089	hKv12.2	2.3
CYL3092	hKv7.4	2.7
CYL3096	hKv7.4/Kv7.5	2.7
CYL3097	hnAChR α 7/ric3	3.0
CYL3099	hKir6.2/SUR2A (KATP)	3.0
CYL3106	hnAChR α 4/ β 2	3.8
CYL3107	hnAChR α 4/ α 6/ β 2	3.8

Document Revision History

Revision Number	Date Released	Revision Details
0	March 2021	<ul style="list-style-type: none">• New document
1.0	July 2023	<ul style="list-style-type: none">• First Revision
2.0	October 2023	<ul style="list-style-type: none">• Second Revision

Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit discoverx.com/legal-policies/ for limited use label license agreement, and trademark information for this product.