

PrecisION[®] hTRPV3 Recombinant Stable Cell Line

Catalog Number CYL3065

Lot Number

See Vial

Contents 2 Vials, 2 x 10⁶ to 4 x 10⁶ in 1 mL

Background Information

The TRPV3 channel belongs to the transient receptor potential channel (TRP) super-family that consists of 7 subfamilies. The channels exist as tetramers where each subunit consists of 6 putative trans-membrane domains. In this respect they have a similar topology to many other voltage-gated channels. The vast majority are permeable to both monovalent cations and calcium (Clapham et al., 2001). The hTRPV3 channel was first cloned in 2002 by two groups (Xu et al., 2002, Smith et al., 2002) based on its homology to known TRPV channels. Additional information can be found on page 2.

Product Information

Description Recombinant HEK 293 cell line expressing the human TRPV3 ligand-gated ion channel (Transient receptor potential cation channel subfamily V member 3, vanilloid receptor-like 3)

Family TRP

Target TRPV3

	Target Protein	Accession Number
1	TRPV3	AJ487035
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

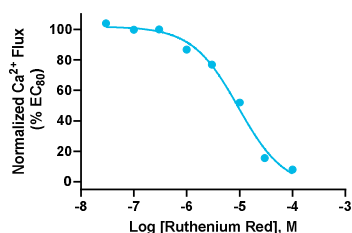
Host Cell Type HEK 293

Application Electrophysiology assay (conventional patch clamp and fluorescent plate-based platforms)

Storage Vials are to be stored in vapor phase of liquid nitrogen

Functional Performance

HEK293 cells expressing hTRPV3 were characterized in terms of their pharmacological and biophysical properties using FLIPR calcium assay.



Electrophysiology Method FLIPR

Reference Agonist 2-APB

Reference Antagonist Ruthenium Red

Antagonist IC₅₀ (μM) 9.42

Passage Stability

This cell line has been confirmed to be stable through at least 12 passages with no significant drop in assay window or change in pharmacology.

Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

Notes

Additional functional (pharmacological and electrophysiological) validation on multiple platforms is available upon request.

Additional Ligand Information

Control Compound Ruthenium Red

Vendor Name : Sigma-Aldrich

Vendor Catalog No. R2751

Additional Background Information

It shares 40-50% homology to TRPV1 (Levine and Alessandri-Haber, 2007) and is activated at temperatures $\geq 34^{\circ}\text{C}$ as well as by 2-Aminoethoxydiphenyl borate (2-APB, Hu et al., 2004) and monoterpenes such as camphor (Vogt-Eisele et al., 2007). In humans it is expressed in skin keratinocytes, trigeminal ganglia, spinal cord and brain (Xu et al., 2002). It is also located in DRG neurons (Xu et al., 2002, Smith et al., 2002, Facer et al., 2007) where it may even form heteromultimers with TRPV1 (Smith et al., 2002). The specific distribution of this channel, coupled with its thermal sensitivity and potential activation by a variety of inflammatory mediators, have suggested that it may represent a novel drug target for the treatment of inflammatory pain.

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com

Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com

General product information: www.discoverx.com

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/license for a list of products that are governed by limited use label license terms and relevant patent and trademark information.