

## PrecisION<sup>®</sup> hTRPV1 Recombinant Stable Cell Line

**Catalog Number** CYL3063

**Lot Number**

See Vial

**Contents** 2 Vials, 2 x 10<sup>6</sup> to 4 x 10<sup>6</sup> in 1 mL

### Background Information

The TRPV1 (Transient Receptor Potential Vanilloid subtype 1) channels cation channels mediate neuronal responses to many noxious stimuli, including heat, low pH, neuropeptides and chemicals e.g. the capsaicin molecule responsible for the sensation of heat from chillies. This receptor may represent a novel target in the treatment of inflammatory pain.

### Product Information

**Description** Recombinant HEK 293 cell line expressing the human TRPV1 ligand-gated ion channel (capsaicin receptor / vanilloid receptor (VR1))

**Family** TRP

**Target** TRPV1

	Target Protein	Accession Number
1	TRPV1	AF196175
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

**Species** Human

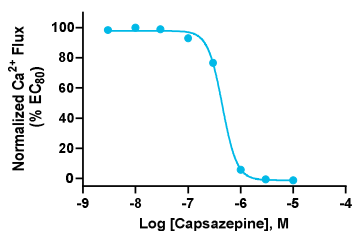
**Host Cell Type** HEK 293

**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 hTRPV1 were characterized in terms of their pharmacological and biophysical properties using FLIPR calcium assay.



**Electrophysiology Method** FLIPR

**Reference Agonist** Capsaicin

**Reference Antagonist** Capsazepine

**Antagonist IC<sub>50</sub> (μM)** 0.45

### 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** Capsazepine

**Vendor Name :** Tocris

**Vendor Catalog No.** 0464

### Additional Background Information

N/A

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)

**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

**General product information:** [www.discoverx.com](http://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](http://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.