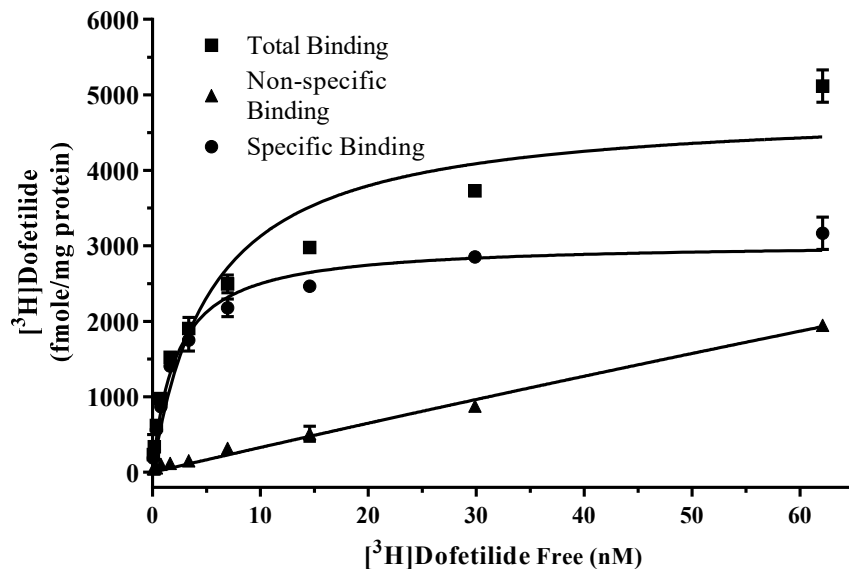


**PRODUCT DATASHEET**
**Precision™ Recombinant hERG Potassium Ion Channel Membrane Preparation**

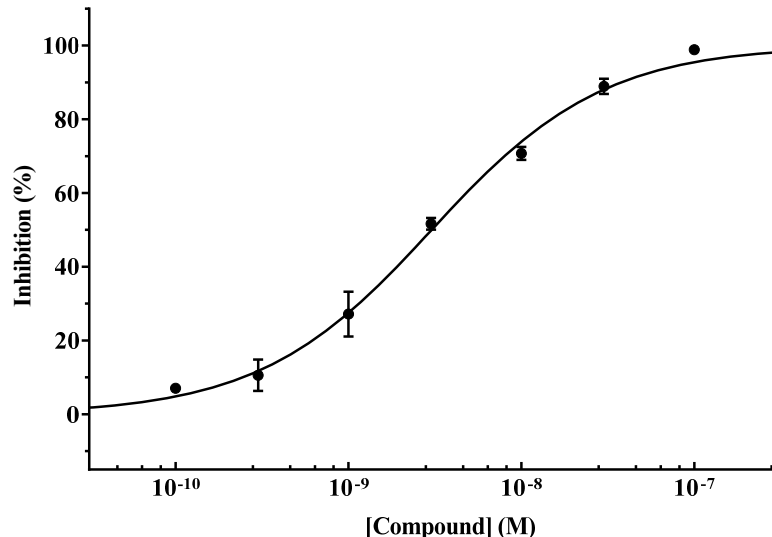
<b>CATALOG NUMBER:</b>	CYL4039	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	22K1204A	<b>VOLUME/CONCENTRATION:</b>	1 mL, 2 mg/mL

**BACKGROUND:** The human ether-a-go-go related gene (hERG) is a potassium ion channel which is essential for normal cardiac repolarization. In drug screening models, the hERG K<sup>+</sup> channel has been indicated to inhibit a wide variety of compounds, and its blockage can lead to cardiac QT interval prolongation and life threatening arrhythmias (Murphy *et al.* 2006). Cardiac safety relating to I<sub>Kr</sub> K<sup>+</sup> channels has become a major concern of regulatory agencies, as hERG channel inhibition has been identified as the firmest link to QT prolongation (Chiu *et al.* 2004). Eurofins hERG membrane preparations are crude membrane preparations made from HEK293 stable recombinant cell lines (Eurofins # CYL3039), which are ideal HTS tools for screening antagonists against the hERG channel. The membrane preparations exhibit a K<sub>d</sub> of 2.5 nM for [<sup>3</sup>H]-Dofetilide. Typically, with 10 μg/well hERG Membrane Prep and 3.0 nM [<sup>3</sup>H]-Dofetilide, a ≥3-fold signal-to-background ratio is obtained.

**APPLICATIONS:** Radioligand binding assay



**Figure 1. Saturation binding for hERG.** 10 μg/well hERG Membrane Preparation was incubated with increasing amount of <sup>3</sup>H-labeled Dofetilide in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled Dofetilide. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for hERG.** hERG Membrane Preparation at 10 µg/well was incubated with 3.0 nM [<sup>3</sup>H]-Dofetilide and increasing concentrations of unlabeled Dofetilide. Typically ≥3-fold signal: background is obtained.

**SPECIFICATIONS:** 1 unit = 10 µg  
 $B_{max}$ : 3.0 pmol/mg  
 $K_d$ : 2.2 nM  
 Signal:background: ≥3-fold

**TRANSFECTION:** Human ERG (Accession number U04270)

**Species:** Human

**HOST CELLS:** HEK-293

**RECOMMENDED ASSAY CONDITIONS:** Membranes are mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a nonbinding 96-well plate, and incubated for 1-2 h. Prior to filtration, an FC 96-well harvest plate (EMD Millipore cat. # MAHF C1H) is coated with 0.33% polyethyleneimine for 30 min. Binding reaction is transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The plate is dried and counted.

**Binding buffer:** 10mM HEPES, pH 7.4, 130mM NaCl, 5mM KCl, 0.8mM MgCl<sub>2</sub>, 1mM NaEGTA, 10mM glucose, 0.1% BSA, filtered and stored at 4°C.

**Radioligand:** [<sup>3</sup>H]-Dofetilide (PerkinElmer, NET-1144)

**Wash Buffer:** 25mM Tris, pH 7.4, 130mM NaCl, 5mM KCl, 0.8mM CaCl<sub>2</sub> 0.1% BSA, filtered and stored at 4°C.

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 3-fold signal: background with <sup>3</sup>H-labeled Astemizole.

**PRESENTATION:**

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.

Packaging method: Membrane proteins were adjusted to the indicated concentration in 1 ml packaging buffer, rapidly frozen, and stored at -80°C.

**STORAGE/HANDLING:** Store at -70°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

**REFERENCES:**

1. Ackerman MJ *et al.* (1998) The long QT syndrome: ion channel diseases of the heart. *Mayo Clin. Proc.* 73: 250-269.
2. Chiu PJS *et al.* (2004) Validation of a [<sup>3</sup>H]-Astemizole Binding Assay in HEK293 Cells Expressing hERG K<sup>+</sup> Channels. *J. Pharmacol. Sci.* 95: 311-319.
3. Chouabe C *et al.* (1998) HERG and KvLQT1/IsK, the cardiac K<sup>+</sup> channels involved in long QT syndromes, are targets for calcium channel blockers. *Mol. Pharmacol.* 54: 695-703.
4. Murphy SM *et al.* (2006) Evaluation of functional and binding assays in cells expressing either recombinant or endogenous hERG channel. *J. Pharmacol. Toxicol. Methods* 54: 42-55.
5. Rampe D *et al.* (1997) A mechanism for the proarrhythmic effects of cisapride (Propulsid): high affinity blockade of the human cardiac potassium channel HERG. *FEBS Lett.* 417: 28-32.

**FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION**

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

**Eurofins Pharma Bioanalytics Services US Inc.** is an independent member of Eurofins Discovery Services