

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT β 1 ADRENERGIC RECEPTOR**

PRODUCT INFORMATION

Catalog Number: DC1437-1

Lot Number: DC1437-1-022114

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human ADRB1 cDNA (GenBank Accession Number AF169007.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: 1-2 days after thawing

Data sheet

Background: Norepinephrine is implicated in a wide range of physiological processes through activation of nine different G-protein-coupled receptors (α 1a, α 1b, α 1d, α 2a, α 2b, α 2c, β 1, β 2, β 3). The human β 1-adrenergic receptor is a 477-amino acid protein found in various heart and brain tissues. β 1 has an important role in the contractile action of valves in cardiac and digestive systems.

Application: Functional assays

Figure 1

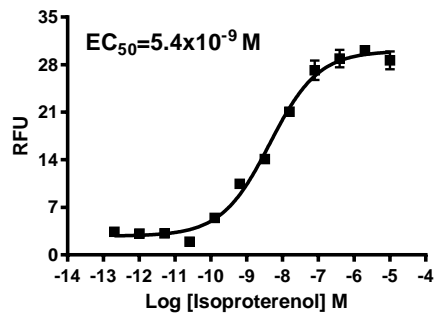


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Oostendorp *et al.* (2000) Contribution of beta-adrenoceptor subtypes to relaxation of colon and oesophagus and pacemaker activity of ureter in wildtype and beta(3)-adrenoceptor knockout mice. *Br J Pharmacol* 130:747-758.

Sato *et al.* (1996) Molecular characterization of pharmacological properties of T-0509 for beta-adrenoceptors. *Eur J Pharmacol* 315:363-367.

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.