

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT β 2 ADRENERGIC RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1438-1a

Lot Number: DC1438-1a-021016

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human ADRB2 cDNA (GenBank Accession Number NM_000024) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 μ g/mL puromycin

Stability: 1-2 days after thawing

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 2-adrenergic receptor was the first 7-transmembrane receptor for a hormone or neurotransmitter to have its crystal structure solved. It has been suggested that the β 2-adrenoceptor may form homodimers as well as oligomers with other receptors. The β 2-adrenoceptor mediates the actions of catecholamines in multiple tissues. They are responsible for relaxation of vascular, uterine, and airway smooth muscle, and are involved in metabolic and endocrine functions.

Application: Functional assays

Figure 1

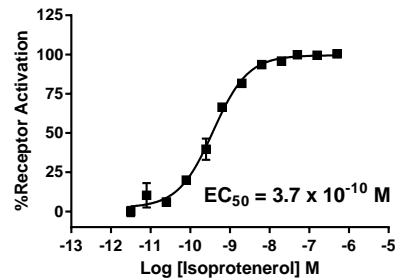


Figure 1. Dose-dependent stimulation of intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Kobilka *et al.* (1987) cDNA for the human beta 2-adrenergic receptor: a protein with multiple membrane-spanning domains and encoded by a gene whose chromosomal location is shared with that of the receptor for platelet-derived growth factor. *Proc Natl Acad Sci USA* 84:46-50.

Frielle *et al.* (1989) Properties of the beta 1- and beta 2-adrenergic receptor subtypes revealed by molecular cloning. *Clin Chem* 35:721-725.

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.