

MULTISCREEN™ β -ARRESTIN2 STABLE CELL LINE HUMAN RECOMBINANT β 2 ADRENERGIC RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CA1438BA2

Lot Number: CA1438BA2-090221

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

Freeze Medium: CellBanker 2 (Amsbio 11891)

Host cell: HEK293T β -Arrestin2

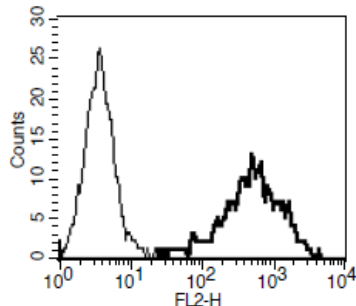
Transfection: Expression vector containing full-length human ADRB2 cDNA (GenBank Accession Number NM_000024) with FLAG tag sequence at N-terminus and ARR2 cDNA (GenBank Accession Number NM_004313.3)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin, 50 μ g/mL hygromycin

Stability: Stable for a minimum of 2 months in continuous culture

Figure 3



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

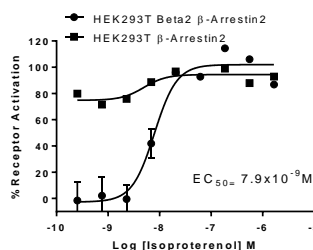


Figure 2

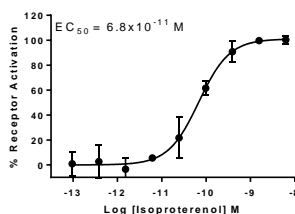


Figure 1. Dose-dependent stimulation from arrestin recruitment upon treatment with ligand, measured with MULTISCREEN™ β -Arrestin Assay Kit (Multispan MSBA01).

Figure 2. Dose-dependent stimulation of intracellular cAMP level upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.