

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT β3 ADRENERGIC RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1439a

Lot Number: DC1439a-050126

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: HEK293T

Transfection: Expression vector containing full-length human ADRB3 cDNA (GenBank Accession Number NM_000025.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Background: The β3-adrenoceptor is also known as the atypical beta-adrenoceptor. It was considered as a potential anti-obesity target by increasing energy expenditure and lipolysis in brown and white adipose tissues of rodents. The functional receptor has been demonstrated in a variety of other tissues including blood vessel, gastrointestinal tract, muscle, and retinal endothelial cells, and its agonist evokes relaxation in some of these tissues such as gall bladder, stomach, small intestine, prostate, colon and bladder.

Application: Functional assays

Figure 1

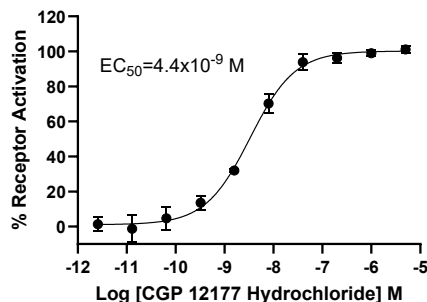


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Emorine *et al.* (1989) Molecular characterization of the human beta 3-adrenergic receptor. *Science* 245:1118-1121.

Fujimura *et al.* (1999) Expression and possible functional role of the beta3-adrenoceptor in human and rat detrusor muscle. *J Urol.* 161:680-685.

FOR RESEARCH USE ONLY.

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