

MULTISCREEN™ β -ARRESTIN2 DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT H4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DCA1030BA2

Lot Number: DCA1030BA2-101221

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

Freeze Medium: CellBanker2

Host cell: HEK293T

Transfection: Expression vector containing full-length human HRH4 cDNA (GenBank accession number NM_021624.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Background: Histamine is one of the most studied biomolecules in medicine and is most notably known for its effects on smooth muscle contraction, vascular permeability and regulation of stomach acid. The histamine receptor H4 has been shown to have a role in chemotaxis and mediator release in a variety of immune cells, such as mast cells, eosinophils, dendritic cells, and T cells. The development of potent H4 receptor antagonists has great potential to open up the pathway for new therapeutic treatments in chronic inflammatory diseases, such as bronchial asthma, allergic gastrointestinal disease, and atopic dermatitis.

Application: Functional assays

Figure 1

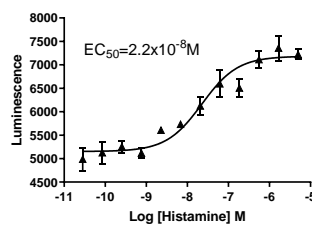


Figure 1. Dose-dependent stimulation from arrestin recruitment upon treatment with ligand, measured with MULTISCREEN™ β -Arrestin Assay Kit (Multispan MSBA01).

References:

Daugherty (2004) Histamine H₄ antagonism: a therapy for chronic allergy?. *Br J Pharmacol* 142:5-7.

Nguyen *et al.* (2001) Discovery of a novel member of the histamine receptor family. *Mol Pharmacol* 59:427-433.

FOR RESEARCH USE ONLY.

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