

MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT H1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: Cm1027

Lot Number: C1027-061714

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length mouse HRH1 cDNA (GenBank Accession Number NM_008285) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/ml puromycin

Stability: Stable in culture for minimum of two months

Background: Histamine H1 receptor (H1 or HRH1) is a subclass of histamine receptors. It mediates the contraction of smooth muscles, increase in capillary permeability due to contraction of terminal venules, and catecholamine release from adrenal medulla, and neurotransmission in the central nervous system. Antihistamines including H1 receptor antagonists are among the most frequently used pharmacologic agents because of a high incident rate of upper respiratory allergies. Clinical and laboratory evidence on histamine and its actions suggests that it has a pathophysiologic role in asthma. These findings have renewed interest in the potential therapeutic role of H1-receptor antagonists, such as, fexofenadine, which has been investigated as potential therapeutic targets for this disease.

Application: Functional assay

Figure 1.

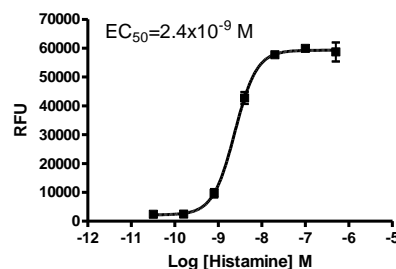


Figure 2.

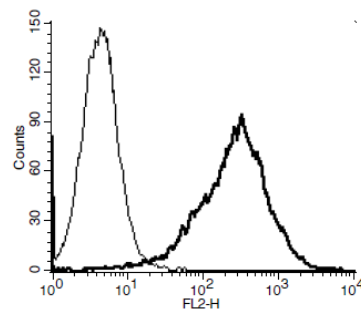


Figure 1. Dose-dependent calcium flux upon treatment with ligand, monitored with FLIPR 384. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Cuss (1999) Beyond the histamine receptor: effect of antihistamines on mast cells. *Clin Exp Allergy* 29 Suppl 3:54-59

Gelfand (2002) Role of histamine in the pathophysiology of asthma: immunomodulatory and anti-inflammatory activities of H1-receptor antagonists. *Am J Med* 113 Suppl 9A:2S-7S.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.