

## MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT H1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** Cm1027

**Lot Number:** C1027-061714

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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

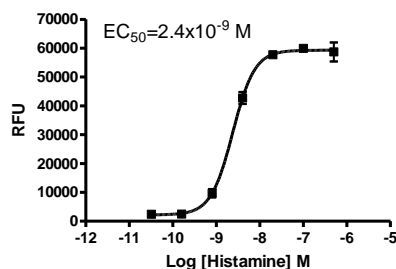
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/ml puromycin

**Stability:** Stable in culture for minimum of two months

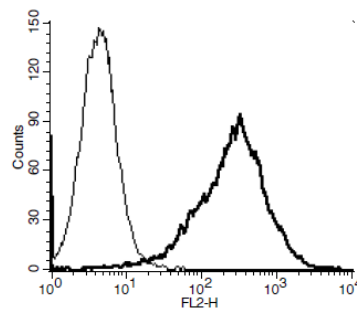
**Background:** Histamine H1 receptor (H1 or HRH1) is a subclass of histamine receptors. It mediates the contraction of smooth muscles, increase in capillary permeability due to contraction of terminal venules, and catecholamine release from adrenal medulla, and neurotransmission in the central nervous 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, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **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.**

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