

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT H2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1028

Lot Number: MC1028-041421

Quantity: 1 vial (10.6 mg/mL),
1mg

Packaging Buffer: 20mM Gly-Gly, 1
mM MgCl₂, 25mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Expression vector
containing full-length human HRH2
cDNA (GenBank accession number
NM_022304.2) with FLAG tag
sequence at N-terminus

Recommended Storage: Liquid
nitrogen upon receiving. Avoid
repeated freeze-thaw

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. H₂ receptors are positively coupled to adenylate cyclase via G_s. It increases the intracellular Ca²⁺ concentrations and release Ca²⁺ from intracellular stores by coupling to G_q. H₂ receptors have been found to be located in a variety of tissues, including the brain, gastric cells, and cardiac tissue. Histamine H₂ receptors have a potent effect on gastric acid secretion, and the inhibition of this secretory process by H₂ receptor antagonists has provided evidence for an important physiological role of histamine in the regulation of gastric secretion. It also regulates gastrointestinal motility and intestinal secretion and is thought to be involved in regulating cell growth and differentiation. It has also been demonstrated to control the relaxation of smooth muscles.

Figure 1

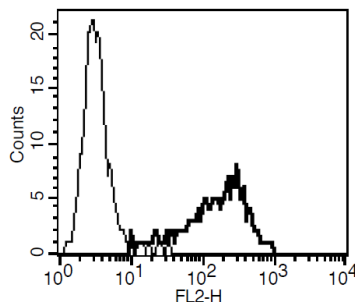


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Hill, S.J. *et al.* (1997) Classification of Histamine Receptors. *Pharmacological Reviews* vol 49 no. 3 253-278

Martínez-Mir, M. I *et al.* (1992) Effect of histamine and histamine analogues on human isolated myometrial strips. *Br J Pharmacol*, 107: 528-531.

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.