

**MULTISCREEN™ MEMBRANES**  
**HUMAN RECOMBINANT H2 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** MC1028

**Lot Number:** MC1028-03292012

**Quantity:** 1 vial (13.1mg/ml)

**Host cell:** HEK293T

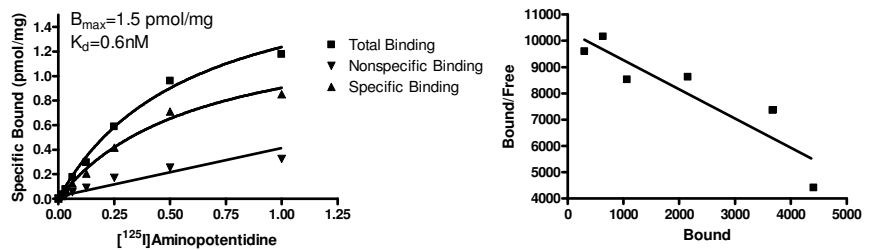
**Transfection:** Expression vector containing full-length human HRH2 cDNA (GenBank accession number NM\_022304.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen

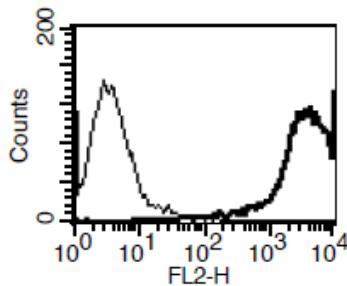
**Background:** Histamine is one of the most studied biomolecules in medicine and is most notably known for its effects on smooth muscle contraction, vascular permeability and regulation of stomach acid. H<sub>2</sub> receptors are positively coupled to adenylate cyclase via G<sub>s</sub>. It increases the intracellular Ca<sup>2+</sup> concentrations and release Ca<sup>2+</sup> from intracellular stores by coupling to G<sub>q</sub>. H<sub>2</sub> receptors have been found to be located in a variety of tissues, including the brain, gastric cells, and cardiac tissue. Histamine H<sub>2</sub> receptors have a potent effect on gastric acid secretion, and the inhibition of this secretory process by H<sub>2</sub> 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.

**Application**  
**Radioligand Binding Assay**

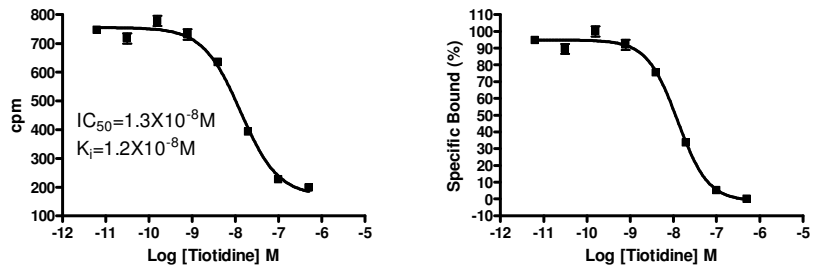
**Figure 1**



**Figure 3**



**Figure 2**



**Figure 1. Saturation binding assay curve:** Specific binding obtained with H<sub>2</sub> membrane with [<sup>125</sup>I]Aminopentidine. Scatchard plot showing one binding site. **Figure 2. Competition binding assay curve:** Dose dependent competition of tiotidine for [<sup>125</sup>I]Aminopentidine 0.05nM. **Figure 3.** 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.**

© 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.

***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](http://www.multispaninc.com)  
[sales@multispaninc.com](mailto:sales@multispaninc.com)  
[support@multispaninc.com](mailto: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.