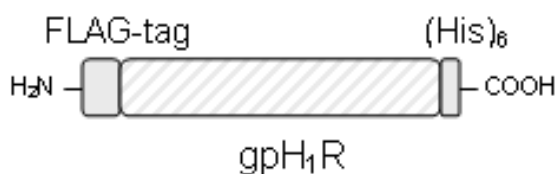


**gpH<sub>1</sub>R + RGS4**

Histamine H<sub>1</sub>-receptor + Regulator of G-protein Signaling 4  
guinea pig, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-1321	1 ml

**For in vitro use only!**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -80 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Molecular Weight:** 60 + 25 kDa

**Accession number:** D15074

**Form:** Membrane suspension (Supplied in 75 mM Tris-HCl pH 7.4, 12.5 mM MgCl<sub>2</sub> and 1 mM EDTA)

**Description:**

The histamine H<sub>1</sub>-receptor is expressed in numerous cell types and couples to G<sub>q</sub>-proteins to activate phospholipase C. In Sf9 insect cells, gpH<sub>1</sub>R couples to the endogenous insect cell G<sub>q</sub>-protein. Co-expression with RGS proteins facilitates detection of high-affinity GTPase activity.

gpH<sub>1</sub>R reconstituted with RGS4 was prepared from Sf9 cells coinfecting with rH<sub>1</sub>R- and RGS4-encoding baculoviruses.

gpH<sub>1</sub>R contains a N-terminal FLAG-tag™ and a C-terminal hexahistidine (His6)-tag for immunological detection, to allow purification and to provide additional protection against proteolysis.

**Selected References:**

Seifert *et al.* (2013) Molecular and cellular analysis of human histamine receptor subtypes. *Trends Pharmacol. Sci.* **34**:33.

Strasser *et al.* (2013) Species-dependent activities of G-protein-coupled receptor ligands: lessons from histamine receptor orthologs. *Trends Pharmacol. Sci.* **34**:13.

Strasser *et al.* (2008) Pharmacological profile of histaprodifens at four recombinant histamine H1 receptor species isoforms. *Pharmacol. Exp. Ther.* **324**:60.