

Product Information



GPR35 Polyclonal Antibody

Catalog No. 10007660

Contents:	This vial contains 50 µg of peptide affinity-purified IgG in 500 µl TBS, pH 7.4, containing 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Synonyms:	G Protein-Coupled Receptor 35, GPCR35
Host:	Rabbit
Antigen:	Synthetic peptide from human GPR35 amino acids 295-304 (AKAHKSQDSL); the antigen alignment with other known sequences is as follows: Human A K A H K S Q D S L Canine A K A H r t Q D S L Murine n t p H K S Q D S q Rat s t p H K S Q D t q
Cross Reactivity:	(+) Human, porcine, and murine GPR35; other species not tested.
Stability:	≥1 year at -20°C
Applications:	The recommended starting dilution for western blot is 1:200 (0.5 µg/ml) and 1:100 (1 µg/ml) for immunohistochemistry (paraffin-embedded sections). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Many G protein-coupled receptor 35 (GPR35) ligands have recently been identified, including kynurenic acid and zaprinast.^{1,2} Multiple isoforms of this receptor have been reported and GPR35 mRNAs were detected in several biological sources including intestine, lymphocytes, skeletal muscle, pancreatic β cells, and tumor cell lines.^{1,3} Cayman's GPR35 polyclonal antibody has been used successfully to detect this receptor in human and murine intestine samples as well as in human, murine, and porcine lymphocytes at 30 kDa on immunoblot.

References

1. Wang, J., Simonavicius, N., Wu, X., *et al.* Kynurenic acid as a ligand for orphan G protein-coupled receptor GPR35. *J. Biol. Chem.* **281(31)**, 22021-22028 (2006).
2. Taniguchi, Y., Tonai-Kachi, H., and Shinjo, K. Zaprinast, a well-known cyclic guanosine monophosphate-specific phosphodiesterase inhibitor, is an agonist for GPR35. *FEBS Lett.* **580**, 5003-5008 (2006).
3. Okumura, S., Baba, H., Kumada, T., *et al.* Cloning of a G-protein-coupled receptor that shows an activity to transform NIH3T3 cells and is expressed in gastric cancer cells. *Cancer Science* **95(2)**, 131-135 (2004).

Related Products

GPR35 Blocking Peptide - Cat. No. 10007661 • 3,4-DAA - Cat. No. 10007980 • Tryptophan Hydroxylase Polyclonal Antibody - Cat. No. 10009396 • Tryptophan Hydroxylase (Phospho-Ser⁵⁸) Polyclonal Antibody - Cat. No. 10009397 • Tryptophan Hydroxylase (Phospho-Ser²⁶⁰) Polyclonal Antibody - Cat. No. 10009398 • Glycine Receptor Polyclonal Antibody - Cat. No. 10009399 • PSD-95 Polyclonal Antibody - Cat. No. 10009506 • NMDA Receptor NR2B Subunit (Phospho-Tyr¹²⁵²) Polyclonal Antibody - Cat. No. 10009759 • NMDA Receptor NR2B Subunit (Phospho-Tyr¹³³⁶) Polyclonal Antibody - Cat. No. 10009760 • NMDA Receptor NR2B Subunit (Phospho-Tyr¹⁴⁷²) Polyclonal Antibody - Cat. No. 1000976

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 01/27/2010



Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com