

PRODUCT INFORMATION



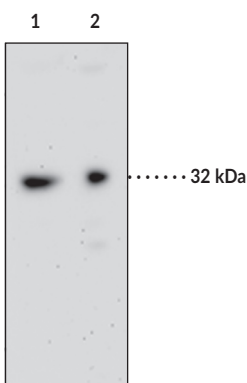
GPR35 Polyclonal Antibody

Item No. 10007660

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	G Protein-Coupled Receptor 35, KYNA Receptor
Immunogen:	Synthetic peptide from the C-terminal region of human GPR35
Species Reactivity:	(+) Human, mouse, and porcine; other species not tested
Uniprot No.:	Q9HC97
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) (paraffin-embedded tissue) and Western blot (WB); the recommended starting dilution for IHC (paraffin-embedded tissue) is 1:100 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Jurkat cell lysate (50 µg)
Lane 2: U937 cell lysate (50 µg)

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/16/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Many G protein-coupled receptor 35 (GPR35) ligands have recently been identified, including kynurenic acid (Item No. 16792) and zaprinast (Item No. 10010421).^{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 mouse intestine samples as well as in human, mouse, 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).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM