

Product Information



Fatty Acid Amide Hydrolase Polyclonal Antibody

Item No. 101600 • Lot. No. XXXXX

- Synonyms:** Anandamide Amidohydrolase; FAAH; Oleamide Hydrolase
- Contents:** This vial contains (100-500 µg of peptide affinity-purified IgG, *lot specific*) in 500 µl TBS, pH *lot specific*, containing *lot specific* mg/ml glycerol, *lot specific*% BSA, and 0.02% sodium azide. This is sufficient for ten (10 x 10 cm) immunoblots.
- Host:** Rabbit
- Antigen:** Synthetic peptide from rat fatty acid amide hydrolase (FAAH) amino acids 561-579 (CLRFMREVEQLMTPQKQPS) conjugated to KLH.¹
Homology of peptide antigen with FAAH from other species:
Rat C L R F M R E V E Q L M T P Q K Q P S
Mouse C L R F M R E V E r L M T P e K r P S
Human C L R F M R E V E r L M T P e K Q s S
Porcine C L R F M R E V E r L M a P g r Q P S
- Cross Reactivity:** (+) Human, rat, mouse, and ferret² FAAH; other species not tested
- Stability:** ≥1 year at -20°C
- Applications:** This polyclonal antibody can be used for detection of FAAH by western blot at a dilution of *lot specific: lot specific*. Immunohistochemistry has not been evaluated and the best working concentration of the antibody should be determined empirically.
- Concentration:** Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 µl final working volume for western blotting.

Laboratory Procedures

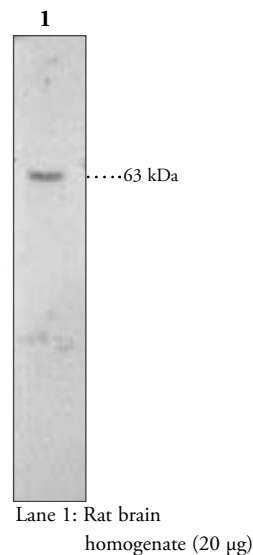
FAAH catalyzes the hydrolysis of biologically significant fatty acid amides.³ Characterization of FAAH, its substrates, and inhibitors have helped to partially elucidate molecular regulation of sleep,⁴ nociception,^{5,6} and cancer.⁷ FAAH is an intracellular enzyme linked to the plasma-membrane *via* its N-terminal domain.^{8,9} The purified enzyme from rat has an estimated molecular mass of 63,000.⁸ Cloning and characterization of porcine, mouse, and human FAAH indicate highly conserved homology with the rat sequence.^{9,10} Northern and immunoblot analyses reveals that FAAH exists in a wide variety of tissues and is particularly abundant in liver, pancreas, brain, testes, uterus, small intestine, and ocular tissue.^{2,3,9,11}

References

1. Tsou, K., Nogueron, M.I., Muthian, S., *et al. Neurosci. Lett.* **254**, 137-140 (1998).
2. Van Sickle, M.D., Oland, L.D., Ho, W., *et al. Gastroenterology* **121**, 767-774 (2001).
3. Fowler, C.J., Jonsson, K.-O., and Tiger, G. *Biochem. Pharmacol.* **62**, 517-526 (2001).
4. Boger, D.L., Henriksen, S.J., and Cravatt, B.F. *Current Pharmaceutical Design* **4**, 303-314 (1998).
5. Cravatt, B.F., Demarest, K., Patricelli, M.P., *et al. Proc. Natl. Acad. Sci. USA* **98(16)**, 9371-9376 (2001).
6. Di Marzo, V., Bisogno, T., Melck, D., *et al. FEBS Lett.* **436**, 449-454 (1998).
7. Bisogno, T., Katayama, K., Melck, D., *et al. Eur. J. Biochem.* **254**, 634-642 (1998).
8. Cravatt, B.F., Giang, D.K., Mayfield, S.P., *et al. Nature* **384**, 83-87 (1996).
9. Giang, D.K. and Cravatt, B.F. *Proc. Natl. Acad. Sci. USA* **94**, 2238-2242 (1997).
10. Goparaju, S.K., Kurahashi, Y., Suzuki, H., *et al. Biochim. Biophys. Acta* **1441**, 77-84 (1999).
11. Katayama, K., Ueda, N., Kurahashi, Y., *et al. Biochim. Biophys. Acta* **1347**, 212-218 (1997).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/101600



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, 12/20/2011