

# Product Information



## Fatty Acid Amide Hydrolase Polyclonal Antibody

Catalog No. 101600 • Lot. No. XXXXX

- Synonyms:** Anandamide Amidohydrolase; FAAH; Oleamide Hydrolase
- Contents:** This vial contains XXX µg peptide affinity-purified IgG in XXX µl TBS, pH XX, containing XX mg/ml glycerol, XX% 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 (CLRFRMREVEQLMTPQKQPS) conjugated to KLH.<sup>1</sup>  
Homology of peptide antigen with FAAH from other species:  
Rat CLRFRMREVEQLMTPQKQPS  
Murine CLRFRMREVE r LMTP e K r P S  
Human CLRFRMREVE r LMTP e K Q s S  
Porcine CLRFRMREVE r L M a P g r Q P S
- Cross-reactivity:** (+) Human, rat, murine, and ferret<sup>2</sup> 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 X:XXX. Immunohistochemistry has not been evaluated and the best working concentration of the antibody should be determined empirically.

### Laboratory Procedures

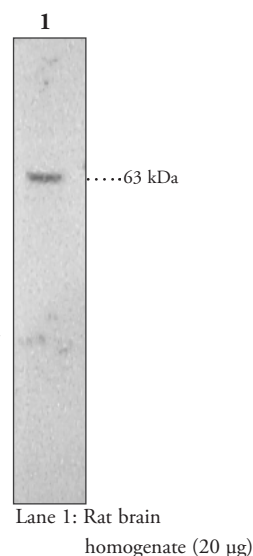
FAAH catalyzes the hydrolysis of biologically significant fatty acid amides.<sup>3</sup> Characterization of FAAH, its substrates, and inhibitors have helped to partially elucidate molecular regulation of sleep,<sup>4</sup> nociception,<sup>5,6</sup> and cancer.<sup>7</sup> FAAH is an intracellular enzyme linked to the plasma-membrane *via* its N-terminal domain.<sup>8,9</sup> The purified enzyme from rat has an estimated molecular mass of 63,000.<sup>8</sup> Cloning and characterization of porcine, murine, and human FAAH indicate highly conserved homology with the rat sequence.<sup>9,10</sup> 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.<sup>2,3,9,11</sup>

### References

1. Tsou, K., Noguero, 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

Methyl Arachidonoyl Fluorophosphonate - Cat. No. 70660 • Arachidonoyl Ethanolamide - Cat. No. 90050 • Decanoyl *p*-Nitroaniline - Cat. No. 90349 • 9-Octadecenamide - Cat. No. 90375 • CB<sub>1</sub> Receptor Polyclonal Antibody - Cat. No. 101500 • CB<sub>2</sub> Receptor Polyclonal Antibody - Cat. No. 101550 • Fatty Acid Amide Hydrolase Blocking Peptide - Cat. No. 301600 • Arachidonoyl Ethanolamide-d<sub>8</sub> - Cat. No. 390050



### 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 NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

#### 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 under separate cover to the MSD supervisor at your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

Copyright Cayman Chemical Company, 12/10/2007