

# Product Information

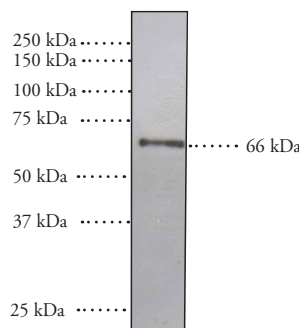


## Fatty Acid Amide Hydrolase Western Ready Control

Catalog No. 10010182

<b>Synonyms:</b>	Anandamide Amidohydrolase, FAAH, Oleamide Hydrolase
<b>Source:</b>	Rat recombinant hexahistidine-tagged protein expressed in <i>E. coli</i>
<b>M<sub>r</sub>:</b>	~66 kDa (as provided), 63 kDa (native)
<b>Purity:</b>	Whole cell lysate; a single band is detected on a western blot when probed with Cayman's Fatty Acid Amide Hydrolase Polyclonal Antibody (Catalog No. 101600)
<b>Supplied in:</b>	Laemmli buffer (2% SDS, 10% glycerol, 62.5 mM Tris-HCl, pH6.8, 5% 2-mercaptoethanol, bromophenol blue)
<b>Stability:</b>	≥2 years at -20°C
<b>Applications:</b>	Positive control for western blotting

Fatty acid amide hydrolase (FAAH) catalyzes the hydrolysis of biologically significant fatty acid amides.<sup>1</sup> Characterization of FAAH, its substrates, and inhibitors have helped to partially elucidate the molecular regulation of sleep,<sup>2</sup> nociception,<sup>3,4</sup> and cancer.<sup>5</sup> FAAH is an intracellular enzyme linked to the plasma-membrane *via* its N-terminal domain.<sup>6,7</sup> The purified enzyme from rat has an estimated molecular mass of 63,000.<sup>6</sup> Cloning and characterization of pig, murine, and human FAAH indicate highly conserved homology with the rat sequence.<sup>7</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>1,7-9</sup>



Western blot of rat recombinant FAAH probed with Cayman's Fatty Acid Amide Hydrolase Polyclonal Antibody (Catalog No. 101600).

### Laboratory Procedures

Western Ready Controls are designed to serve as positive controls on western blots. This FAAH Western Ready Control is supplied pre-boiled in 55 µl of Laemmli buffer. To use, thaw and load 10 µl of FAAH Western Ready Control per well and follow a standard western blotting protocol. The amount needed per well will depend on the antibody and detection system used. Due to the presence of the 6X His-tag on the FAAH provided in this control, the protein will migrate as a band ~3 kDa higher than the native, non-tagged protein.

### References

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

### Related Products

Fatty Acid Amide Hydrolase Polyclonal Antibody - Cat. No. 101600 • Fatty Acid Amide Hydrolase Blocking Peptide - Cat. No. 301600

**WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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 under separate cover to the MSDS supervisor at 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, 06/20/2008

### 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