

Product Information

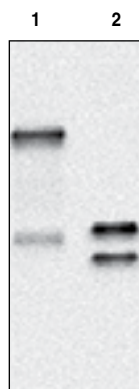


ANGPTL3 (mouse) Monoclonal Antibody (Clone Kairos3-3741)

Item No. 10807

Contents:	This vial contains a 1 mg/ml solution of antibody in PBS, pH 7.4.
Synonyms:	ANG-5, Angiotensin-5, Angiotensin-Like 3, Angiotensin-Related Protein 3, ANGPT5, FHBL2
Antigen:	Recombinant mouse ANGPTL3
Host:	Rat, clone Kairos3-3741
Isotype:	IgG _{2ak}
Cross Reactivity:	(+) Mouse ANGPTL3; (-) mouse ANGPTL4 and human ANGPTL3; rat and other species not tested
Stability:	≥6 months at -20°C
Applications:	ELISA and western blot (WB); the recommended starting dilution for ELISA (direct or indirect) is 1:2,000-1:5,000 and 1:1,000-1:5,000 for WB. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

ANGPTL3 regulates angiogenesis and also directly regulates lipid, glucose, and energy metabolism. ANGPTL3 (mouse) monoclonal antibody (clone Kairos3-3741) recognizes mouse ANGPTL3. It detects a band of ~69 kDa by WB and cross-reacts with mouse ANGPTL3 fibrinogen-like domain (FLD).



Lane 1: Recombinant ANGPTL3 (mouse)
Lane 2: Recombinant ANGPTL3 (mouse) (FLD)

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10807

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, 08/24/2012

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