

Product Information

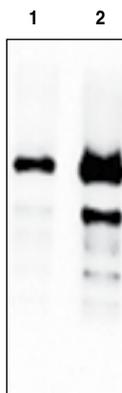


FTO (mouse) Monoclonal Antibody (Clone FT 342-1)

Item No. 10815

Contents:	This vial contains a 1 mg/ml solution of antibody in PBS, pH 7.4.
Synonyms:	Fat Mass- and Obesity-associated Protein, FATSO
Antigen:	Recombinant mouse FTO
Host:	Rat, clone FT 342-1
Isotype:	IgG _{2ak}
Cross Reactivity:	(+) Mouse and rat FTO; (-) human FTO.
Stability:	≥6 months at -20°C
Applications:	ELISA and western blot (WB). The recommended starting dilution for ELISA (direct or indirect) and WB: 1:2,000-1:5,000.

Fat mass- and obesity-associated gene (FTO) is the gene responsible for mouse 'fused toes' mutation. An association between FTO genotype and type 2 diabetes has been confirmed. The presence of the FTO rs9939609A-allele was found to be positively correlated with other symptoms of the metabolic syndrome, including higher fasting insulin, glucose, and triglycerides, and lower HDL-cholesterol.



Lane 1: Recombinant rat FTO
Lane 2: Recombinant murine FTO

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10815

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, 12/14/2011