

Product Information



Nampt Monoclonal Antibody (Clone OMNI 379)

Item No. 10813

| | |
|--------------------------|---|
| Contents: | This vial contains a 1 mg/ml solution of antibody in PBS, pH 7.4. |
| Synonyms: | Nicotinamide Phosphoribosyltransferase, PBEF1, Pre-B Cell Colony Enhancing Factor 1, Visfatin |
| Antigen: | Recombinant human Nampt |
| Host: | Mouse, clone OMNI 379 |
| Isotype: | IgG _{2ak} |
| Cross Reactivity: | (+) Human, mouse, and rat Nampt |
| Stability: | ≥6 months at -20°C |
| Applications: | ELISA, flow cytometry (FC), immunocytochemistry (ICC), immunohistochemistry (IHC), immunoprecipitation (IP), and western blot (WB). The recommended starting dilution for ELISA (direct or indirect) is 1:2,000-1:10,000; 1:1,000 for FC, ICC, IHC, and IP; and 1:2,000-1:5,000 for WB. |

Nampt is a 52 kDa adipokine secreted by adipose tissue that is involved in the biosynthesis of nicotinamide adenine dinucleotide (NAD⁺). Two forms of Nampt exist, an intracellular form (iNampt) and an extracellular form (eNampt). While the function of iNampt as an essential and rate-limiting NAD⁺ biosynthetic enzyme is well established, the physiological role of eNampt is still a matter of debate. Nampt has various functions, including the promotion of vascular smooth muscle cell maturation and inhibition of neutrophil apoptosis. It activates the insulin receptor and has insulin-mimetic effects, lowering blood glucose and improving insulin sensitivity. The protein is highly expressed in visceral fat and serum levels of the protein correlate with obesity.



Lane 1: Recombinant human Nampt (His-tagged)
Lane 2: Recombinant human Nampt (FLAG[®]-tagged)
Lane 3: Recombinant mouse Nampt (His-tagged)
Lane 4: Recombinant rat Nampt (His-tagged)

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10813

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.

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 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, 08/26/2013