

CERTIFICATE of ANALYSIS



Nampt Monoclonal Antibody (Clone OMNI 379)

Item No. 10813 • Batch No. 0810417 • Expires 24MAY2028

Overview

Formulation:	A 1 mg/ml solution in PBS, pH 7.4
Concentration:	(no concentration)
Storage/stability:	-20°C/1yr
Applications:	ELISA, FC, ICC, IHC, IP, and WB
Species reactivity:	(+)Human Nampt(+)Mouse Nampt(+)Rat Nampt

Reviewed and approved by: Emily Tucker

Approved date: 04AUG2025

A handwritten signature in black ink, appearing to read "Emily Tucker".

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM