

CERTIFICATE of ANALYSIS



PTP1B (human, recombinant)

Item No. 10010896 • Batch No. 0803486 • Expires 02APR2027

Overview

Formulation:	100 mM Tris, pH 7.2, with 100 mM sodium chloride, 1 mM EDTA, and 25% glycerol
Concentration:	0.67 mg/mL
Storage/stability:	-80°C/1yr
Specific activity:	18.00 U/mg
Unit definition:	One unit of enzyme produces 1 μ mol of p-nitrophenol/min at 25°C in 100 mM sodium citrate, pH 6.0, containing 10 mM p-nitrophenyl phosphate.
Activity:	12.00 U/ml
Purity:	≥90% estimated by SDS-PAGE

Reviewed and approved by: Beth Hurst

Approved date: 03APR2026

Elizabeth Hurst

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