

CERTIFICATE of ANALYSIS



SARS-CoV-2 nsp13 Helicase (E. coli expressed)

Item No. 30589 • Batch No. 0676378 • Expires 30JAN2025

Overview

Formulation:	50 mM Tris-HCl, pH 7.5, with 5 μ M zinc chloride, 1 mM DTT, 1 mM EDTA, and 15% glycerol
Concentration:	0.56 mg/mL
Storage/Stability:	-80°C/1yr
Specific activity:	485,378.00 U/mg
Unit definition:	One unit is defined as the amount of enzyme required to release 1 pmol of fluorescein-labeled ssDNA from a fork DNA substrate containing a quencher on the complimentary strand per minute at 25°C under the following conditions: 20 mM HEPES, pH 7.5, 20 mM NaCl, 5% glycerol, 5 mM MgCl ₂ , 0.01% BSA, 2 mM DTT, 2 μ M unlabeled complementary ssDNA, 1 mM ATP, and 200 nM forked DNA substrate.
Activity:	271,811.70 U/ml
Purity:	≥70% estimated by SDS-PAGE

Reviewed and approved by: Carly Norris

Approved date: 13APR2023

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