

PRODUCT INFORMATION



Histone H3 Peptide Substrate (1-21)

Item No. 10877

Overview

Contents: A lyophilized peptide
Synonym: H3 Peptide
Peptide Sequence: ARTKQTARKSTGGKAPRKQLA
FW: 2,254.6

Properties

Form: Solid
Storage: -20°C (as supplied)
Stability: ≥3 years
Purity: ≥95% by HPLC

Application

Water or aqueous buffers: At least 20 mg/ml

Other methods were not tested, therefore optimal working concentration/dilution should be determined empirically.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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, 11/08/2023

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