

# 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:** ≥1 year  
**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, 09/10/2015

#### 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