

PRODUCT INFORMATION



Casein Kinase 2 Substrate (trifluoroacetate salt)

Item No. 24848

Formal Name:	L-arginyl-L-arginyl-L-arginyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-threonyl-L- α -glutamyl-L- α -glutamyl-L-glutamic acid, trifluoroacetate salt	
Synonym:	CK2 Substrate	
MF:	$C_{52}H_{87}N_{19}O_{24} \cdot XCF_3COOH$	H-Arg-Arg-Arg-Glu-Glu-Glu-Thr-Glu-Glu-Glu-OH
FW:	1,362.4	$\cdot XCF_3COOH$
Purity:	$\geq 95\%$	
Supplied as:	A lyophilized powder	
Storage:	$-20^\circ C$	
Stability:	≥ 4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Casein kinase 2 substrate (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the casein kinase 2 substrate (trifluoroacetate salt) in water. The solubility of casein kinase 2 substrate (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Casein kinase 2 substrate is a peptide substrate for casein kinase 2 (CK2).¹ It is selectively phosphorylated by CK2 over a panel of eight kinases and the amount of phosphorylated peptide can be used to quantify CK2 enzyme activity.

Reference

1. Kuenzel, E.A. and Krebs, E.G. A synthetic peptide substrate specific for casein kinase II. *Proc. Natl. Acad. Sci. U.S.A.* **82(3)**, 737-741 (1985).

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, 11/18/2022

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