

PRODUCT INFORMATION



Malantide (trifluoroacetate salt)

Item No. 40150

Formal Name:	L-arginyl-L-threonyl-L-lysyl-L-arginyl-L-serylglycyl-L-seryl-L-valyl-L-tyrosyl-L- α -glutamyl-L-prolyl-L-leucyl-L-lysyl-L-isoleucine, trifluoroacetate salt	
Peptide Sequence:	RTKRSGSVYEPLKI-OH	H-Arg-Thr-Lys-Arg-Ser-Gly-Ser-Val-Tyr-Glu-Pro-Leu-Lys-Ile-OH
MF:	C ₇₂ H ₁₂₄ N ₂₂ O ₂₁ • XCF ₃ COOH	
FW:	1,633.9	
Purity:	≥98%	• XCF ₃ COOH
Ex./Em. Max:	285/302 nm	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

Malantide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the malantide (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Malantide is a synthetic peptide derived from the β -subunit of human phosphorylase kinase and is a fluorescent substrate for PKA and PKC.^{1,2} Upon phosphorylation by PKA or PKC, the intrinsic tyrosine fluorescence decreases and can be used to quantify PKA and PKC activity. Malantide displays an emission at 302 nm upon excitation at 285 nm.

References

1. Malencik, D.A. and Anderson, S.R. Characterization of a fluorescent substrate for the adenosine 3',5'-cyclic monophosphate-dependent protein kinase. *Anal. Biochem.* **132(1)**, 34-40 (1983).
2. Zhao, Z.H., Malencik, D.A., and Anderson, S.R. Characterization of a new substrate for protein kinase C: Assay by continuous fluorometric monitoring and high performance liquid chromatography. *Biochem. Biophys. Res. Commun.* **176(3)**, 1454-1461 (1991).

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, 02/20/2024

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