

PRODUCT INFORMATION



LRRKtide (trifluoroacetate salt)

Item No. 37461

Formal Name:	L-arginyl-L-leucylglycyl-L-arginyl-L- α -aspartyl-L-lysyl-L-tyrosyl-L-lysyl-L-threonyl-L-leucyl-L-arginyl-L-glutamyl-L-isoleucyl-L-arginyl-L-glutamine, trifluoroacetate salt	
Synonyms:	Membrane-organizing Extension Spike Protein (550-564), Moesin (550-564)	H-Arg-Leu-Gly-Arg-Asp-Lys-Tyr-Lys-Thr-Leu-Arg-Gln-Ile-Arg-Gln-OH • XCF ₃ COOH
Peptide Sequence:	RLGRDKYKTLRQIRQ-OH	
MF:	C ₈₃ H ₁₄₇ N ₃₁ O ₂₂ • XCF ₃ COOH	
FW:	1,931.3	
Purity:	≥98%	
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

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

Description

LRRKtide is a peptide substrate for leucine-rich repeat kinase 2 (LRRK2), an enzyme frequently mutated in patients with Parkinson's disease, that corresponds to amino acids 550-564 of membrane-organizing extension spike protein (moesin).¹ It has been used to identify inhibitors of LRRK2.

Reference

1. Leveridge, M., Collier, L., Edge, C., *et al.* A high-throughput screen to identify LRRK2 kinase inhibitors for the treatment of Parkinson's Disease using RapidFire mass spectrometry. *J. Biomol. Screen.* **21(2)**, 145-155 (2016).

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, 10/13/2023

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