

PRODUCT INFORMATION

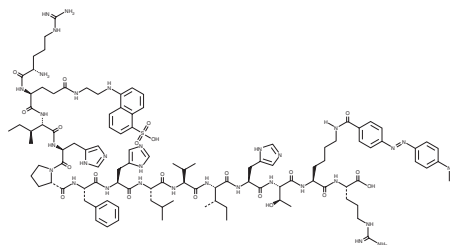


Renin Fluorogenic Substrate

Item No. 10006906

Overview and Properties

Contents:	This vial contains 1 mg of a crystalline solid.
CAS Registry No.:	791068-69-4
Formal Name:	L-arginyl-N-[2-[(5-sulfo-1-naphthalenyl)amino]ethyl]-L-glutamyl-L-isoleucyl-L-histidyl-L-prolyl-L-phenylalanyl-L-histidyl-L-leucyl-L-valyl-L-isoleucyl-L-histidyl-L-threonyl-N ⁶ -[4-[[4-(dimethylamino)phenyl]azo]benzoyl]-L-lysyl-L-arginine
Uniprot No.:	P00797
MF:	C ₁₀₉ H ₁₅₆ N ₃₂ O ₂₁ S
FW:	2,282.7
Form:	Solid
Purity:	≥95%
Ex./Em. Max:	340/490 nm
Storage:	-20°C (as supplied)
Stability:	≥3 years



Laboratory Procedures

The renin fluorogenic substrate consists of the normal peptide substrate for renin which has been linked to the fluorophore EDANS at one end and to a non-fluorescent quenching molecule (Dabcyl) at the other. After cleavage by renin, the product (peptide-EDANS) is brightly fluorescent and can be easily analyzed using an excitation wavelength of 340 nm and emission wavelengths of 485-510 nm.

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