

PRODUCT INFORMATION



Amyloid- β (1-8) Peptide

Item No. 10241

Overview and Properties

Contents: This vial contains 1 mg of lyophilized peptide.
Synonym: A β (1-8)
Peptide Sequence: Human β -amyloid amino acids 1-8; DAEFRHDS
FW: 975.96
Storage: -20°C (as supplied); use reconstituted aliquots within 30 days if -20°C storage
Stability: \geq 3 years
Solubility: 10 mg/ml with water or aqueous buffers

Description

Amyloid- β (1-8) is a wild-type control for the mutation-containing amyloid- β (1-8, A2V) peptide (Item No. 10229).

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, 11/03/2023

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