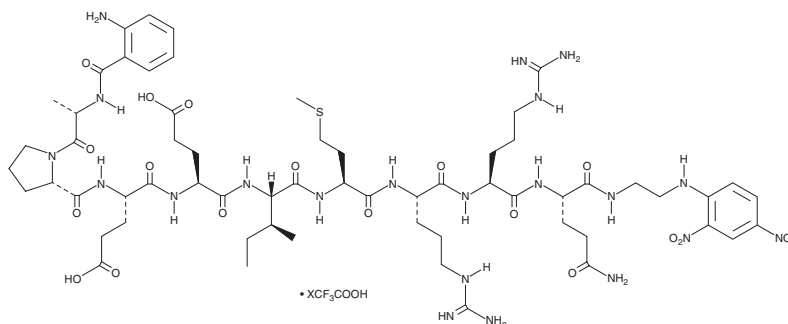


PRODUCT INFORMATION



Abz-Ala-Pro-Glu-Glu-Ile-Met-Arg-Arg-Gln-EDDnp (trifluoroacetate salt) Item No. 16467

Formal Name: N-(2-aminobenzoyl)-L-alanyl-L-prolyl-L- α -glutamyl-L- α -glutamyl-L-isoleucyl-L-methionyl-L-arginyl-L-arginyl-N¹-[2-[(2,4-dinitrophenyl)amino]ethyl]-L-glutamamide, trifluoroacetate salt
MF: C₆₁H₉₃N₂₁O₁₉S • XCF₃COOH
FW: 1,456.6
Purity: ≥95%
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

Abz-Ala-Pro-Glu-Glu-Ile-Met-Arg-Arg-Gln-EDDnp (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Abz-Ala-Pro-Glu-Glu-Ile-Met-Arg-Arg-Gln-EDDnp (trifluoroacetate salt) in water.

Description

Abz-Ala-Pro-Glu-Glu-Ile-Met-Arg-Arg-Gln-EDDnp is a fluorescence-quenched peptide substrate for human neutrophil elastase ($k_{cat}/K_m = 531 \text{ mM}^{-1}\text{s}^{-1}$).¹ Enzymatic peptide hydrolysis disrupts the Abz:EDDnp donor-acceptor pair allowing fluorescence measurement of Abz (ex = 320 nm; em = 420 nm).

Reference

1. Korkmaz, B., Attucci, S., Juliano, M.A., *et al.* Measuring elastase, proteinase 3 and cathepsin G activities at the surface of human neutrophils with fluorescence resonance energy transfer substrates. *Nat. Protoc.* **3**(6), 991-1000 (2008).

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, 11/07/2022

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