

PRODUCT INFORMATION



Bz-IEGR-pNA (acetate)

Item No. 27567

CAS Registry No.: 1325307-57-0

Formal Name: N-benzoyl-L-isoleucyl-L- α -glutamylglycyl-N-(4-nitrophenyl)-L-argininamide, monoacetate

Synonyms: Bz-Ile-Glu-Gly-Arg-pNA,
Bz-Ile-Glu-Gly-Arg-p-nitroanilide,
Factor Xa Chromogenic Substrate

MF: $C_{32}H_{43}N_9O_9 \cdot C_2H_4O_2$

FW: 757.8

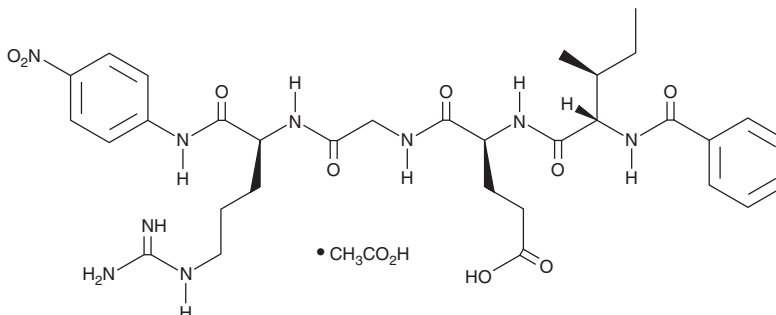
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 224, 314 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years



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

Description

Bz-IEGR-pNA is a colorimetric substrate for Factor Xa.¹ Factor Xa preferentially binds to and cleaves the Ile-Glu-Gly-Arg (IEGR) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of Factor Xa activity.

Reference

1. Aruelli, L., Friberger, P., Karlsson, G., *et al.* A new sensitive and highly specific chromogenic peptide substrate for factor Xa. *Thromb. Res.* **11**(5), 595-609 (1977).

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, 11/02/2022

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