

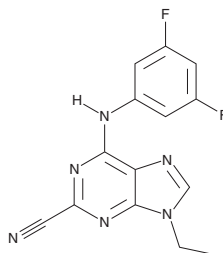
PRODUCT INFORMATION



Cruzain-IN-1

Item No. 37350

CAS Registry No.: 1199523-24-4
Formal Name: 6-[(3,5-difluorophenyl)amino]-9-ethyl-9H-purine-2-carbonitrile
Synonym: Cruzain Inhibitor 1
MF: C₁₄H₁₀F₂N₆
FW: 300.3
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

Cruzain-IN-1 is supplied as a solid. A stock solution may be made by dissolving the cruzain-IN-1 in the solvent of choice, which should be purged with an inert gas. Cruzain-IN-1 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

Cruzain-IN-1 is an inhibitor of the *Trypanosoma* protease cruzain (IC₅₀ = 0.2 nM).¹ It is selective for cruzain over the *Trypanosoma* proteases cathepsin L-like protein and cathepsin B-like protease (IC₅₀s = <6 and 6,000 nM, respectively) and a panel of 71 mammalian cysteine proteases, aspartyl proteases, peptidases, matrix metalloproteinases (MMPs), and serine proteases at 1 μM. Cruzain-IN-1 is active against *T. brucei* (MIC₅₀ = 25 μM).

Reference

1. Mott, B.T., Ferreira, R.S., Simeonov, A., *et al.* Identification and optimization of inhibitors of Trypanosomal cysteine proteases: Cruzain, rhodesain, and TbCatB. *J. Med. Chem.* **53**(1), 52-60 (2010).

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, 04/03/2024

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