

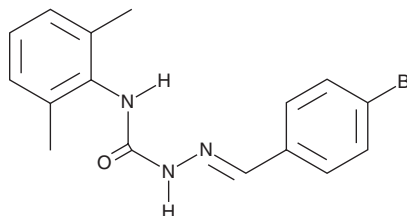
PRODUCT INFORMATION



EGA

Item No. 21882

CAS Registry No.: 415687-81-9
Formal Name: 2-[(4-bromophenyl)methylene]-
N-(2,6-dimethylphenyl)-
hydrazinecarboxamide
MF: C₁₆H₁₆BrN₃O
FW: 346.2
Purity: ≥98%
UV/Vis.: λ_{max}: 291 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.

Laboratory Procedures

EGA is supplied as a crystalline solid. A stock solution may be made by dissolving the EGA in the solvent of choice, which should be purged with an inert gas. EGA is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

EGA is an inhibitor of endosomal trafficking.¹ It increases accumulation of fluorescently labeled EGF in early endosomes in HeLa cells when used at a concentration of 20 μM. EGA inhibits cell death induced by anthrax lethal toxin in RAW264.7 cells (IC₅₀ = 1.7 μM) as well as by additional acid-dependent bacterial toxins, including diphtheria toxin, *P. aeruginosa* ExoA, and *H. ducreyi* cytolethal distending toxin (Hd-CDT), in various cell types. It also inhibits infection of Vero and HeLa cells by the low pH-dependent lymphocytic choriomeningitis (LCMV; Armstrong strain) and influenza A/WSN/33 viruses, respectively.

Reference

1. Gillespie, E.J., Ho, C.L., Balaji, K., *et al.* Selective inhibitor of endosomal trafficking pathways exploited by multiple toxins and viruses. *Proc. Natl. Acad. Sci. U.S.A.* **110(50)**, E4904-E4912 (2013).

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, 12/02/2022

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