

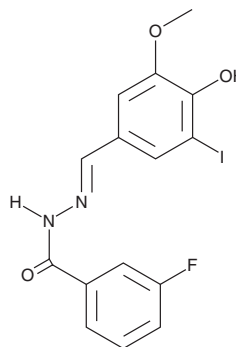
PRODUCT INFORMATION



Endosidin 2

Item No. 21888

CAS Registry No.: 1839524-44-5
Formal Name: 3-fluoro-benzoic acid, (2E)-2-[(4-hydroxy-3-iodo-5-methoxyphenyl)methylene]hydrazide
Synonym: ES2
MF: C₁₅H₁₂FIN₂O₃
FW: 414.2
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 328 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Endosidin 2 (ES2) is supplied as a crystalline solid. A stock solution may be made by dissolving the ES2 in the solvent of choice. ES2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ES2 in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

ES2 is a cell-permeable benzylidene-benzohydrazide that binds to the exocyst component of the 70 kDa (EXO70) subunit of the exocyst complex ($K_d = 253 \mu\text{M}$, EXO70A1).¹ ES2 binding inhibits exocytosis and endosomal recycling in plant and mammalian cells. ES2 disrupts protein trafficking between the endosome and plasma membrane, which enhances trafficking to the vacuole for degradation. It also inhibits recycling of endocytosed transferrin to the plasma membrane in HeLa cells and can target multiple isoforms of mammalian EXO70, resulting in misregulation of exocytosis.

Reference

1. Zhang, C., Brown, M.Q., van de Ven, W., *et al.* Endosidin2 targets conserved exocyst complex subunit EXO70 to inhibit exocytosis. *Proc. Natl. Acad. Sci. USA* **113**(1), E41-E50 (2016).

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, 08/14/2017

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