

PRODUCT INFORMATION



Ferristatin II

Item No. 36621

CAS Registry No.: 1937-37-7

Formal Name: 4-amino-3-[2-[4'-[2-(2,4-diaminophenyl) diazenyl][1,1'-biphenyl]-4-yl]diazenyl]-5-hydroxy-6-(2-phenyldiazenyl)-2,7-naphthalenedisulfonic acid, disodium salt

Synonyms: Chlorazol Black, C.I. 30235, NSC 8679

MF: C₃₄H₂₅N₉O₇S₂ • 2Na

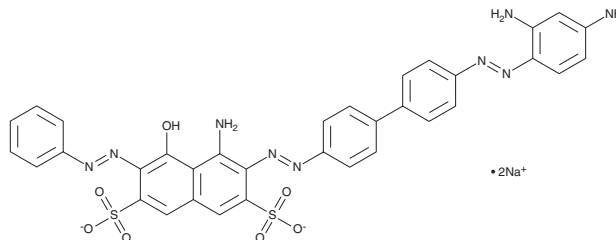
FW: 781.7

UV/Vis.: λ_{max}: 596 nm

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

Ferristatin II is supplied as a solid. Aqueous solutions of ferristatin II can be prepared by directly dissolving the solid in aqueous buffers. The solubility of ferristatin II in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ferristatin II is an inhibitor of iron uptake.¹ It inhibits iron uptake (IC₅₀ = ~12 μM), as well as induces degradation of transferrin receptor protein 1 (TfR1), an effect that can be reversed by the lipid membrane disruptor nystatin, in HeLa cells. *In vivo* ferristatin II (0.2, 10, and 40 mg/kg) reduces serum iron levels and transferrin saturation in rats. It also suppresses traumatic brain injury-induced lipid peroxidation, neuronal apoptosis, and the number of cortical iron deposits in rats.²

References

1. Byrne, S.L., Buckett, P.D., Kim, J., *et al.* Ferristatin II promotes degradation of transferrin receptor-1 *in vitro* and *in vivo*. *PLoS One* **8**(7), e70199 (2013).
2. Cheng, Y., Qu, W., Li, J., *et al.* Ferristatin II, an iron uptake inhibitor, exerts neuroprotection against traumatic brain injury via suppressing ferroptosis. *ACS Chem. Neurosci.* **13**(5), 664-675 (2022).

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, 04/17/2025

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