

PRODUCT INFORMATION



TC HSD 21

Item No. 29730

CAS Registry No.: 330203-01-5

Formal Name: 5-[(3-bromo-4-hydroxyphenyl)methylene]-3-(4-methoxyphenyl)-2-thioxo-4-thiazolidinone

Synonym: TC Hydroxysteroid Dehydrogenase 21

MF: $C_{17}H_{12}BrNO_3S_2$

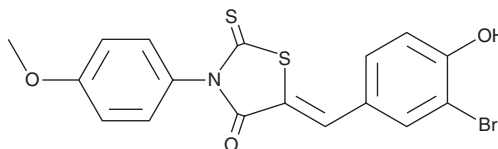
FW: 422.3

Purity: $\geq 98\%$

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

TC HSD 21 is supplied as a solid. A stock solution may be made by dissolving the TC HSD 21 in the solvent of choice, which should be purged with an inert gas. TC HSD 21 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

TC HSD 21 is an inhibitor of 17β -hydroxysteroid dehydrogenase type 3 (17β -HSD3; $IC_{50} = 14$ nM in HeLa cells expressing the recombinant human enzyme).¹ It is selective for 17β -HSD3 over 17β -HSD1, 17β -HSD2, estrogen receptor α (ER α), the androgen receptor, and the glucocorticoid receptor.

Reference

1. Harada, K., Kubo, H., Tanaka, A., *et al.* Identification of oxazolidinediones and thiazolidinediones as potent 17β -hydroxysteroid dehydrogenase type 3 inhibitors. *Bioorg. Med. Chem. Lett.* **22**(1), 504-507 (2012).

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, 10/21/2022

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