

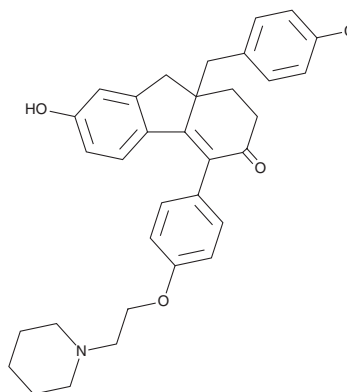
PRODUCT INFORMATION



CMP8

Item No. 36585

CAS Registry No.: 851107-28-3
Formal Name: 9a-[(4-chlorophenyl)methyl]-1,2,9,9a-tetrahydro-7-hydroxy-4-[4-[2-(1-piperidinyl)ethoxy]phenyl]-3H-fluoren-3-one
MF: C₃₃H₃₄ClNO₃
FW: 528.1
Purity: ≥98%
UV/Vis.: λ_{max}: 346 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

CMP8 is supplied as a solid. A stock solution may be made by dissolving the CMP8 in the solvent of choice, which should be purged with an inert gas. CMP8 is soluble in DMSO.

Description

CMP8 is a mutant estrogen receptor α (ER α) antagonist and ligand used for bump and hole, a technique to study single protein isoforms when familial homology would interfere with other isoform-selective approaches.¹ It is selective for an ER α triple mutant containing a lysine-to-methionine substitution at position 384 (ER α ^{L384M}), a methionine-to-glycine substitution at position 421 (ER α ^{M421G}), and a glycine-to-arginine substitution at position 521 (ER α ^{G521R}), over wild-type ER α and ER β (IC₅₀s = 29, 1,085, and 2,232 nM, respectively). It reduces reporter gene expression in HeLa cells expressing the triple mutant ER α (EC₅₀ = 41 nM).

Reference

1. Gallinari, P., Lahm, A., Koch, U., et al. A functionally orthogonal estrogen receptor-based transcription switch specifically induced by a nonsteroid synthetic ligand. *Chem. Biol.* **12**(8), 883-893 (2005).

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, 07/21/2023

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