

PRODUCT INFORMATION

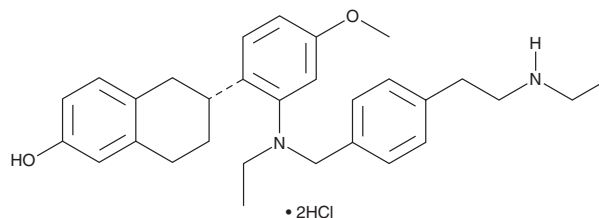


Elacestrant (hydrochloride)

Item No. 37298

CAS Registry No.: 1349723-93-8
Formal Name: 6R-[2-[ethyl[[4-[2-(ethylamino)ethyl]phenyl]methyl]amino]-4-methoxyphenyl]-5,6,7,8-tetrahydro-2-naphthalenol, dihydrochloride

Synonym: RAD901
MF: C₃₀H₃₈N₂O₂ • 2HCl
FW: 531.6
Purity: ≥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

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

Description

Elacestrant is a selective estrogen receptor degrader (SERD).¹ It selectively binds to estrogen receptor α (ER α) over ER β (IC₅₀s = 48 and 870 nM, respectively) and reduces ER α levels in MCF-7 cells (EC₅₀ = 0.6 nM). Elacestrant reduces MCF-7 proliferation induced by 17 β -estradiol (E₂; Item No. 10006315) (EC₅₀ = 4 pM). *In vivo*, elacestrant (30 and 60 mg/kg) reduces E₂-induced tumor growth in an MCF-7 mouse xenograft model. It increases survival in an MCF-7 intracranial mouse xenograft model. Elacestrant (0.1-3 mg/kg) protects against bone loss in ovariectomized rats.

Reference

1. Garner, F., Shomali, M., Paquin, D., *et al.* RAD1901: A novel, orally bioavailable selective estrogen receptor degrader that demonstrates antitumor activity in breast cancer xenograft models. *Anticancer Drugs* 26(9), 948-956 (2015).

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, 09/26/2023

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