PRODUCT INFORMATION



Nilutamide-d₆

Item No. 28694

CAS Registry No.:	1189477-66-4	р Н
Formal Name:	5,5-bis(methyl-d ₃)-3-[4-nitro-3-(trifluoromethyl)	
	phenyl]-2,4-imidazolidinedione	
MF:	$C_{12}H_4D_6F_3N_3O_4$	
FW:	323.3	
Chemical Purity:	≥98% (Nilutamide)	D 0.
Deuterium		
Incorporation:	≥99% deuterated forms (d ₁ -d ₆); ≤1% d ₀	$\langle \rangle$
Supplied as:	A solid	$\rightarrow = \langle$
Storage:	-20°C	E-C NO
Stability:	≥4 years	130 1102

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

Laboratory Procedures

Nilutamide-d₆ is intended for use as an internal standard for the quantification of nilutamide (Item No. 23953) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Nilutamide-d₆ is supplied as a solid. A stock solution may be made by dissolving the nilutamide-d₆ in the solvent of choice, which should be purged with an inert gas. Nilutamide- d_{4} is soluble in DMSO.

Description

Nilutamide is a non-steroidal antiandrogen that is an antagonist of the androgen receptor.¹ It reverses testosterone-induced increases in cell proliferation with an IC50 value of 412 nM in mouse mammary carcinoma cells in vitro.² It has antiandrogenic activity in both intact and castrated rats, inhibiting testosterone propionate-induced prostate growth when administered at a dose of 2.5 mg/kg.³ Formulations containing nilutamide have been used in combination with surgical castration in the treatment of metastatic prostate cancer.

References

- 1. Signh, S.M., Gauthier, S., and Labrie, F. Androgen receptor antagonists (antiandrogens): Structure-activity relationships. Curr. Med. Chem. 7(2), 211-247 (2000).
- 2. Simard, J., Singh, S.M., and Labrie, F. Comparison of in vitro effects of the pure antiandrogens OH-flutamide, Casodex, and nilutamide on androgen-sensitive parameters. Urology 49(4), 580-586 (1997).
- 3. Wakeling, A.E., Furr, B.J.A., Glen, A.T., et al. Receptor binding and biological activity of steroidal and nonsteroidal antiandrogens. J. Steroid Biochem. 15, 355-359 (1981).

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

uyer 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, 12/05/2022

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