

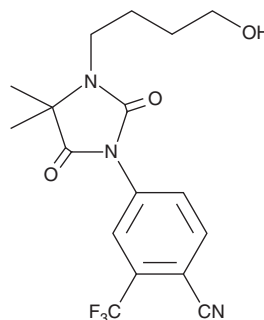
PRODUCT INFORMATION



RU-58841

Item No. 24680

CAS Registry No.: 154992-24-2
Formal Name: 4-[3-(4-hydroxybutyl)-4,4-dimethyl-2,5-dioxo-1-imidazolidiny]-2-(trifluoromethyl)-benzonitrile
MF: C₁₇H₁₈F₃N₃O₃
FW: 369.3
Purity: ≥98%
UV/Vis.: λ_{max}: 262 nm
Supplied as: A crystalline 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

RU-58841 is supplied as a crystalline solid. A stock solution may be made by dissolving the RU-58841 in the solvent of choice, which should be purged with an inert gas. RU-58841 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RU-58841 is approximately 33 mg/ml in ethanol and DMF and approximately 25 mg/ml in DMSO.

RU-58841 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RU-58841 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. RU-58841 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RU-58841 is a non-steroidal antiandrogen that binds to the androgen receptor ($K_d = 1.8, 1.1, 5.6, 1.3,$ and 1.4 nM for human, rat, mouse, hamster prostate, and hamster flank organ receptors, respectively).¹ *In vivo*, topical administration of RU-58841 (1-100 µg per animal) reduces size of the flank organ without affecting sex organs in a dose-dependent manner in hamsters. It has no effect on sex organ weight in intact rats when administered percutaneously, subcutaneously, or orally at doses up to 1 mg per animal. RU-58841 (5% v/v) also induces hair regrowth in the bald frontal scalps of postpubertal stump-tailed macaques, a model for male pattern baldness.²

References

1. Battmann, T., Bonfils, A., Branche, C., *et al.* RU 58841, a new specific topical antiandrogen: A candidate of choice for the treatment of acne, androgenetic alopecia and hirsutism. *J. Steroid Biochem. Mol. Biol.* **48(1)**, 55-60 (1994).
2. Obana, N., Chang, C., and Uno, H. Inhibition of hair growth by testosterone in the presence of dermal papilla cells from the frontal bald scalp of the postpubertal stump-tailed macaque. *Endocrinology* **138(1)**, 356-361 (1997).

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, 12/06/2022

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