

PRODUCT INFORMATION



MK-0773

Item No. 43202

CAS Registry No.: 606101-58-0
Formal Name: (4aS,4bS,6aS,7S,9aS,9bS,11aR)-3-fluoro-2,4a,4b,5,6,6a,7,8,9,9a,9b,10,11,11a-tetradecahydro-N-(3H-imidazo[4,5-b]pyridin-2-ylmethyl)-1,4a,6a-trimethyl-2-oxo-1H-indeno[5,4-f]quinoline-7-carboxamide

Synonym: PF-05314882

MF: C₂₇H₃₄FN₅O₂

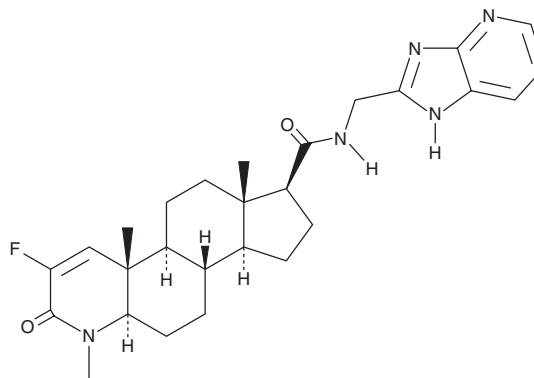
FW: 479.6

Purity: ≥90%

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

MK-0773 is supplied as a solid. A stock solution may be made by dissolving the MK-0773 in the solvent of choice, which should be purged with an inert gas. MK-0773 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

MK-0773 is a selective androgen receptor modulator (SARM).¹ It selectively binds to the androgen receptor (AR; IC₅₀ = 6.6 nM) over the glucocorticoid and progesterone receptors (IC₅₀s = >2 μM for both). It is a partial agonist of AR that increases AR-dependent transcription in a reporter assay using MDA-MB-453 cells endogenously expressing AR. MK-0773 (6 mg/kg) increases the bone formation rate with less than a 5% increase in uterine weight in ovariectomized rats and increases seminal vesicle weight to a lesser extent than dihydrotestosterone (DHT) in orchidectomized rats when administered at a dose of 45 mg/kg. It reduces serum total cholesterol levels but also reduces serum HDL levels in ovariectomized rats when administered at a dose of 25 mg/kg.

Reference

- Schmidt, A., Kimmel, D.B., Bai, C., *et al.* Discovery of the selective androgen receptor modulator MK-0773 using a rational development strategy based on differential transcriptional requirements for androgenic anabolism versus reproductive physiology. *J. Biol. Chem.* **285**(22), 17054-17064 (2010).

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, 04/10/2026

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