

Bicalutamide

Androgen receptor inhibitor

Nonsteroidal antiandrogen. Effective in the treatment of prostate cancer. Induces cell death by a pathway that is independent of changes in mitochondrial membrane potential and Bcl-2 action.

Citations: 22

[View Online »](#)

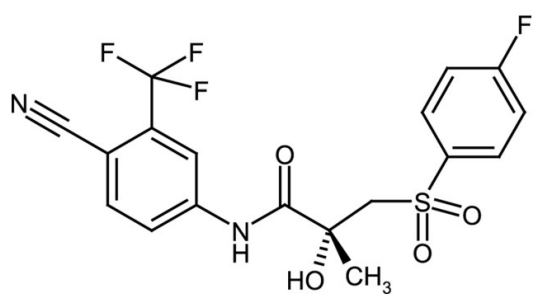
Ordering Information

[Order Online »](#)

ALX-270-476-M100	100mg
ALX-270-476-M500	500mg
ALX-270-476-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(4-Cyano-3-(trifluoromethyl)phenyl)-3-((4-fluorophenyl)sulfonyl)-2-hydroxy-2-methylpropanamide

Appearance White crystalline powder.

CAS 90357-06-5

Couple Target Androgen receptor

Couple Type Inhibitor

Formula $C_{18}H_{14}F_4N_2O_4S$

MW 430.4

Purity ≥98% (HPLC)

Solubility Soluble in acetone and tetrahydrofuran; slightly soluble in 100% ethanol or methanol; practically insoluble in water.

Source Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com