

Fostriecin

Antitumor antibiotic

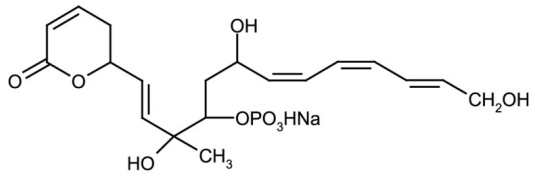
Antitumor antibiotic. Catalytic inhibitor of topoisomerase II. Strong protein phosphatase 2A (PP2A) inhibitor ($IC_{50}=3.2nM$). Weak inhibitor for PP1 ($IC_{50}=131\mu M$). No apparent effect on PP2B. The binding site for fostriecin on PP2A is different from that of okadaic acid (see Prod. No. ALX-350-003).

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year when stored at -20°C. Use fresh solution.
Handling	Protect from light. Packaged under inert gas. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Phosphotrienin, CI-920
Appearance	Colorless solid.
CAS	87860-39-7
Couple Target	Serine/threonine-protein phosphatase, Topoisomerase
Couple Type	Inhibitor
Formula	$C_{19}H_{26}O_9PNa$
MW	452.4
Purity	≥98% (HPLC)
RTECS	UQ0600000
Solubility	Soluble in water (50mg/ml), methanol (10mg/ml) or 100% ethanol (10mg/ml).
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-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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