

DPN

Estrogen receptor β activator

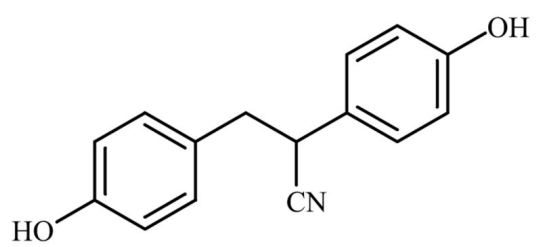
Selective estrogen receptor ER β agonist displaying 70-fold greater selectivity over ER α (EC_{50} = 0.85 and 66nM, respectively). It displays neuroprotective activity in various *in vitro* and *in vivo* models.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Diarylpropionitrile, 2,3-bis(4-Hydroxyphenyl)-propionitrile
Appearance	White solid.
CAS	1428-67-7
Couple Target	Estrogen receptor
Couple Type	Activator, Ligand
Formula	$C_{15}H_{13}NO_2$
MW	239.3
MeltingPoint	197-199°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol (1.5mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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