

AAL-993

VEGFR inhibitor

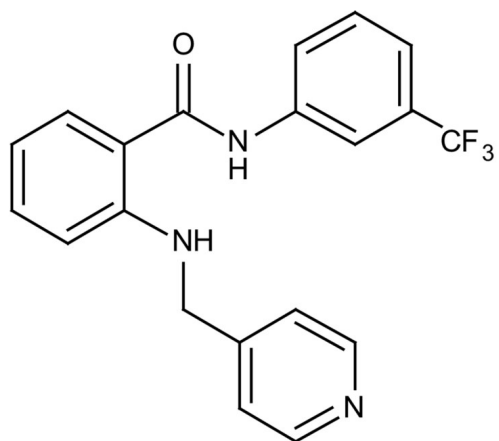
Highly potent inhibitor of VEGFR-1 (IC_{50} =130nM), VEGFR-2 (IC_{50} =23nM) and VEGFR-3 (IC_{50} =18nM). At higher concentrations it inhibits PDGFR- β (640nM), c-Kit (236nM) and CSF-1R (380nM).

Citations: 2

[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 room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2-[(4-Pyridyl)methyl]amino-N-[3-(trifluoromethyl)phenyl]benzamide
Appearance	Pale yellow solid.
CAS	269390-77-4
Couple Target	VEGFR
Couple Type	Inhibitor
Formula	$C_{20}H_{16}F_3N_3O$
MW	371.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25 mg/ml) or ethanol (15 mg/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)