

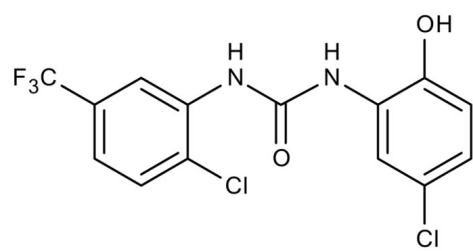
NS 1738

Selective positive allosteric modulator of $\alpha 7$ nAChR.

A selective positive allosteric modulator selective for $\alpha 7$ nicotinic acetylcholine receptors ($\alpha 7$ nAChR). NS 1738 exhibits cognitive enhancing effects in vivo.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	1-(5-chloro-2-hydroxyphenyl)-3-(2-chloro-5-(trifluoromethyl)phenyl)urea, $\alpha 7$ nAChR positive allosteric modulator
Appearance	White crystalline solid.
CAS	501684-93-1
Couple Target	Acetylcholine receptor
Couple Type	Activator
Formula	$C_{14}H_9Cl_2F_3N_2O_2$
MW	365.1
Purity	$\geq 99\%$ (HPLC)
Solubility	Soluble in DMSO or ethanol.

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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