

ORG-12962

Serotonin receptor activator

Selective serotonin 5-HT_{2C} agonist. Displays antiaversive effects in a panic-like anxiety rat model. May display appetite regulating properties similar to other 5-HT_{2C} agonists.

Citations: 3

[View Online »](#)

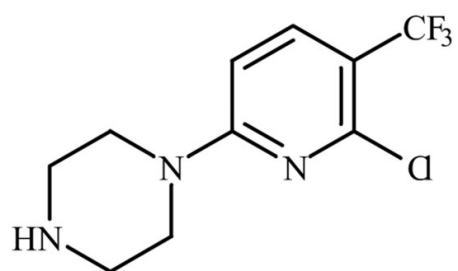
Ordering Information

[Order Online »](#)

BML-NS705-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 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	1-[6-Chloro-5-trifluoromethyl)-2-pyridinyl]-piperazine
Appearance	Pale yellow solid.
CAS	210821-63-9
Couple Target	Serotonin receptor
Couple Type	Activator, Ligand
Formula	$C_{10}H_{11}ClF_3N_3$
Identity	Determined by 1H -NMR.
MW	265.7
Purity	≥95% (NMR)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/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-be@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-
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)