

NS-1643

K⁺ channel activator

A novel and selective HERG K⁺ channel activator (EC₅₀=10.5 µM). Activates ERG2 channels. Displays antiarrhythmic properties in rabbit models.

Citations: 3

[View Online »](#)

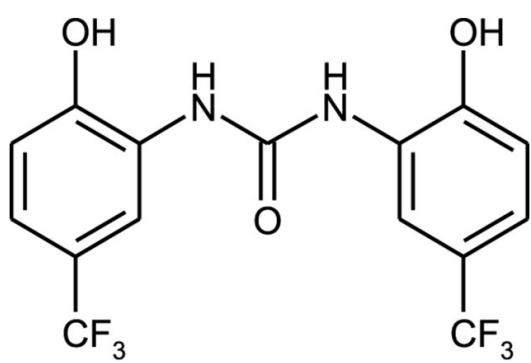
Ordering Information

[Order Online »](#)

BML-KC169-0010	10mg
BML-KC169-0050	50mg

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 after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N,N'-Bis[2-hydroxy-5-(trifluoromethyl)phenyl] urea
Appearance	White solid
CAS	448895-37-2
Couple Target	Potassium channel
Couple Type	Activator
Formula	$C_{15}H_{10}F_6N_2O_3$
MW	380.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 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-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-
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)