

BML-283

Pyruvate kinase M2 activator

Potent ($AC_{50} = 0.111\mu M$) and selective (over PKM1, PKR and PKL) activator of the tumor cell specific M2 isoform of pyruvate kinase (PKM2). Pyruvate kinase is a potential target for cancer chemotherapy. A useful new tool for altering aberrant glucose metabolism in cancer cells (Warburg effect).

Citations: 1

[View Online »](#)

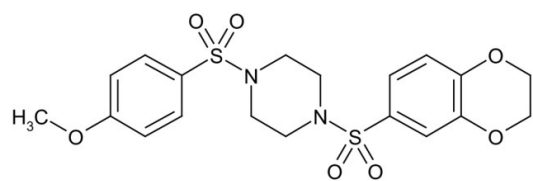
Ordering Information

[Order Online »](#)

BML-EI404-0025	25mg
----------------	------

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 one 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	NCGC00030335, 1-(2,3-Dihydrobenzo[b][1,4]dioxin-6-ylsulfonyl)-4-(4-methoxyphenylsulfonyl)-piperazine
Appearance	White solid.
Couple Target	Pyruvate kinase
Couple Type	Activator
Formula	$C_{19}H_{22}N_2O_7S_2$
MW	454.5
MeltingPoint	229-231°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (12mg/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)