

# STO-609

## CaM kinase kinase inhibitor

STO-609 is a specific CaM kinase kinase (CaM-KK) inhibitor.  $K_i=0.2\mu\text{M}$  and  $0.04\mu\text{M}$  for  $\alpha$  and  $\beta$  isotypes respectively. Demonstrates >80-fold selectivity for CaM-KK over CaMK1, CaMK2, CaMK4, MLCK, PKA, PKC, and p42 MAPK. Concentrations up to  $10\mu\text{g/ml}$  do not affect cellular viability. It is a tool for exploring the role of the CaM-KK pathway in various physiological processes.

Citations: 4

[View Online »](#)

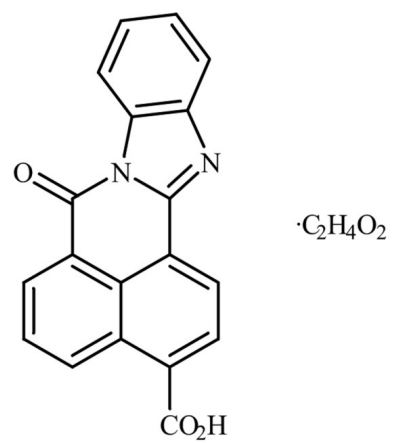
## Ordering Information

[Order Online »](#)

BML-EI389-0005	5mg
BML-EI389-0025	25mg

## 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	7-Oxo-7H-benzimidazo[2,1-a]benz[de]isoquinoline-3-carboxylic acid, acetate salt
Appearance	Yellow solid.
CAS	1173022-21-3
Couple Target	CaM-KK
Couple Type	Inhibitor
Formula	$C_{19}H_{10}N_2O_3 \cdot C_2H_4O_2$
MW	374.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/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)