

# ESI-09

## Membrane-permeant inhibitor of Epac 1 and Epac 2

ESI-09 is a membrane-permeant inhibitor of both exchange proteins activated directly by cyclic AMP, Epac 1 and Epac 2. ESI-09 displays *in vitro* values of 3.2  $\mu\text{M}$  for Epac 1 and 1.4  $\mu\text{M}$  for Epac 2 with 25  $\mu\text{M}$  cAMP as competing activator (Almahariq et al. 2013). It displays an at least 100-fold selectivity for Epac proteins compared to cyclic AMP-dependent protein kinase I and II in assays (Almahariq et al. 2013).

*BLG-B133-25 (5 x 5  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

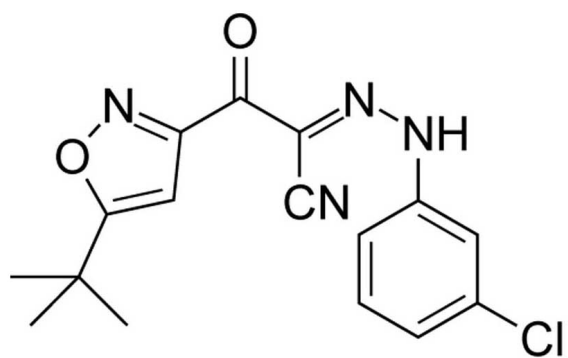
### Ordering Information

[Order Online »](#)

<b>BLG-B133-05</b>	5 $\mu\text{mol}$
<b>BLG-B133-25</b>	5x5 $\mu\text{mol}$

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 3-[5-(tert.-Butyl)isoxazol-3-yl]-2-[2-(3-chlorophenyl)hydrazono]-3-oxopropanenitrile

**CAS** 263707-16-0

**Couple Type** Modified nucleotides

**Formula**  $C_{16}H_{15}ClN_4O_2$

**MW** 330.8

**Purity** > 95% HPLC

**Quantity** 5 µmol = ~1.7 mg

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)