

# IDE-2

## Induces stem cell differentiation

Cell-permeable molecule capable of directing differentiation of mouse and human embryonic stem cells into endodermal cells ( $EC_{50}=223$  nM). Developmental potential of these chemically derived endodermal cells has been shown via continued differentiation into pancreatic progenitor cells and incorporation into developing gut tubes of live mouse embryos. Nodal/Smad signaling is activated indicating IDE-2 functions via TGF- $\beta$ -signaling pathway.

Citations: 1

[View Online »](#)

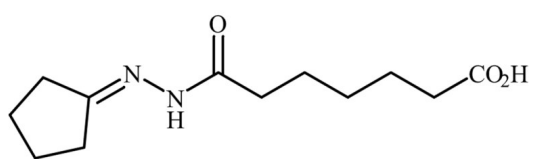
## Ordering Information

[Order Online »](#)

|                |     |
|----------------|-----|
| BML-SC101-0001 | 1mg |
|----------------|-----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                          |  |
|--------------------------|--|
| <b>Use/Stability</b>     | As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to one year. Store solutions at -20°C for up to 3 months. |
| <b>Long Term Storage</b> | -20°C  |
| <b>Shipping</b>          | Blue Ice   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                         |   |
|-------------------------|---|
| <b>Alternative Name</b> | 7-(2-Cyclopentylidenehydrazinyl)-7-oxoheptanoic acid    |
| <b>Appearance</b>       | White solid.  |
| <b>CAS</b>              | 1136466-93-7  |
| <b>Formula</b>          | $C_{12}H_{20}N_2O_3$                                    |
| <b>MW</b>               | 240.3   |
| <b>Purity</b>           | ≥98% (TLC)  |
| <b>Solubility</b>       | Soluble in DMSO (>25 mg/ml) or 100% ethanol (25 mg/ml). |



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)