

# IDE1

## Small molecule inducer of definitive endoderm

IDE1 is a small-molecule compound that promotes the differentiation of embryonic stem cells (ESCs) and induced pluripotent stem cells (iPSCs) into definitive endoderm (DE) lineages. It functions by activating the TGF- $\beta$  signaling pathway, leading to Smad2 phosphorylation and increased Nodal expression. IDE1 has an EC<sub>50</sub> of approximately 125 nM for inducing DE differentiation.

Key features and applications include:

- **Definitive Endoderm Induction:** Efficiently induces Sox17<sup>+</sup>/FoxA2<sup>+</sup> endodermal progenitors from both mouse and human ESCs.
- **TGF- $\beta$  Pathway Activation:** Mimics Activin A by triggering Smad2 phosphorylation, a key step in endoderm specification.
- **Cell-Permeable & Cost-Effective:** Offers a small-molecule alternative to expensive growth factors like Activin A.
- **Versatile in Culture Systems:** Effective in both 2D monolayer and 3D embryoid body (EB) differentiation protocols.
- **Pancreatic Lineage Differentiation:** When combined with FGF10, retinoic acid, or hedgehog inhibitors, IDE1-derived DE cells can further differentiate into Pdx1<sup>+</sup> pancreatic progenitors.

Research Applications:

- Stem cell differentiation protocols
- Developmental biology studies
- Pancreatic and hepatic lineage specification
- High-throughput screening for endoderm inducers

Relevant disease states include:

- **Diabetes Mellitus:** IDE1 is used in protocols to generate pancreatic  $\beta$ -cell precursors, aiding in diabetes research and regenerative therapy development.

- **Liver Disease:** Supports the generation of hepatocyte-like cells from pluripotent stem cells for liver disease modeling and drug screening.
- **Congenital Gut Disorders:** Enables modeling of early gut tube development and related congenital abnormalities.
- **Stem Cell Therapy Research:** Facilitates the development of cell-based therapies by providing a scalable method for generating endodermal derivatives.

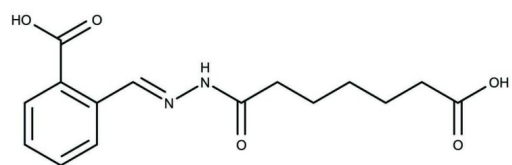
### Ordering Information

[Order Online »](#)

ENZ-CHM352-0025	25mg
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-[(E)-(6-carboxyhexanoylhydrazinylidene)methyl]benzoic acid
Appearance	White solid.
CAS	1160927-48-9
Couple Target	TGF-beta
Couple Type	Activator
Formula	$C_{15}H_{18}N_2O_5$
Identity	Determined by NMR.
MW	306.32
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (≥ 25 mg/mL).

Last modified: July 28, 2025



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)