

6-Chloro-2,3,4,9-tetrahydro-1H-carbazole-1-carboxamide

SIRT inhibitor

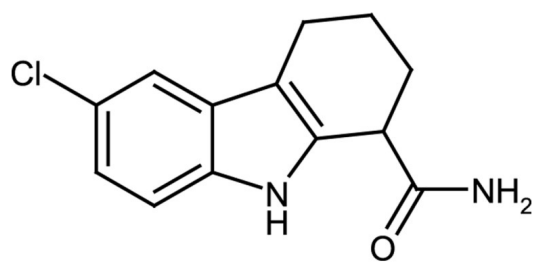
Potent cell permeable and metabolically stable specific inhibitor of hSIRT1 (IC_{50} =98nM *in vivo* / IC_{50} =38nM *in vitro*; compared to hSIRT2: IC_{50} =19 μ M and hSIRT3: IC_{50} =48 μ M) with no effect on human histone deacetylases (HDACs) class I and class II, nor NAD glycohydrolase (IC_{50} >100 μ g). Inhibits the deacetylation of p53 (IC_{50} =1 μ M).

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|---|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 6 months when stored at -20°C in the dark. |
| Handling | Protect from light. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------------|--------------------------------|
| Alternative Name | EX-527 |
| Appearance | White to light yellow powder. |
| CAS | 49843-98-3 |
| Couple Target | SIRT |
| Couple Type | Inhibitor |
| Formula | $C_{13}H_{13}ClN_2O$ |
| Identity | Identity determined by 1H-NMR. |
| MW | 248.7 |
| Purity | ≥95% (HPLC) |
| Purity Detail | Racemic mixture |
| 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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)