

# Caveolin-1 scaffolding domain peptide

Cell permeable eNOS blocker

Cell permeable peptide that blocks eNOS (NOS III) activity and cellular nitric oxide (NO) release *in vitro* and reduces inflammation and tumorigenesis *in vivo*. Caveolin-1 interacts with several lipid-modified signalling ligands, such as EGFR, eNOS, G-protein  $\alpha$ -subunits, PKC $\alpha$ , H-Ras, and Src, via the C1-SD82-101 sequence.

Citations: 9

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-153-064-M001 | 1mg |
| ALX-153-064-M005 | 5mg |

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                                      |
|-------------------|--------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C. |
| Handling          | Protect from light. Hygroscopic.                                                                                                     |
| Long Term Storage | -20°C                                                                                                                                |
| Shipping          | Ambient Temperature                                                                                                                  |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                                                                                                      |
|------------------|------------------------------------------------------------------------------------------------------------------------------------------------------|
| Alternative Name | Cavtratin, AP-Cav, Pen-C1-SD                                                                                                                         |
| Appearance       | White lyophilized powder.                                                                                                                            |
| Formula          | $C_{228}H_{335}N_{61}O_{49}S$                                                                                                                        |
| MW               | 4746.6                                                                                                                                               |
| Purity           | ≥97% (HPLC)                                                                                                                                          |
| Sequence         | H-Arg-Gln-Ile-Lys-Ile-Trp-Phe-Gln-Asn-Arg-Arg-Met-Lys-Trp-Lys-Lys-Asp-Gly-Ile-Trp-Lys-Ala-Ser-Phe-Thr-Thr-Phe-Thr-Val-Thr-Lys-Tyr-Trp-Phe-Tyr-Arg-OH |
| Solubility       | Soluble in DMSO or acetonitrile:water (50:50).                                                                                                       |

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)