

Myelin oligodendrocyte glycoprotein (mouse, rat) (35-55)

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|------------------------------------------------------------------------------------------|
| Alternative Name | MOG (35-55) |
| Appearance | White to off-white powder. |
| CAS | 163913-87-9 |
| Formula | $C_{118}H_{177}N_{35}O_{29}S$ |
| MW | 2582.0 |
| Purity | ≥85% (HPLC) |
| Sequence | H-Met-Glu-Val-Gly-Trp-Tyr-Arg-Ser-Pro-Phe-Ser-Arg-Val-Val-His-Leu-Tyr-Arg-Asn-Gly-Lys-OH |
| Solubility | <1mg/ml soluble in water, methanol, DMSO or dimethyl formamide. |
| Source | Synthetic peptide. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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