

Myelin Oligodendrocyte Glycoprotein (35-55) (rat)

Oligodendrocyte antigen used to induce

5AE
Ordering Information

[Order Online »](#)

LKT-M9645-M001	1mg
LKT-M9645-M002	2mg
LKT-M9645-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MOG (35-55)

CAS 163913-87-9

Formula $C_{118}H_{177}N_{35}O_{29}S$

MW 2581.99

Purity ≥98%

Solubility Soluble in water.



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