

Estradiol Glow, L pack

Fluorescently labeled Estradiol

Ordering Information

[Order Online »](#)

JBS-PR-958L	5x100µg
-------------	---------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance brownish-red, glassy to waxy appearance

Emission Maximum 618 nm

Excitation Maximum 467 nm

Formula $C_{47}H_{67}BrN_2O_2$

Formulation lyophilised

MW 771.95

Purity > 97 % (DC, NMR)

Solubility High: MeOH, EtOH, PrOH, DMSO, DMF Moderate: acetone, CH₂Cl₂, CH₃Cl, THF Low: H₂O and aqueous buffer systems (after presolution in DMSO >1 µM) Insoluble: Et₂O, hydrocarbons such as hexane, heptane, benzene

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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