

Rosiglitazone

Thiazolidinedione; PPAR γ agonist.

Rosiglitazone stimulates AMP-activated protein kinase.

Ordering Information

[Order Online »](#)

LKT-R5773-M500	500mg
LKT-R5773-G001	1g
LKT-R5773-G005	5g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BRL-49653

CAS 122320-73-4

Couple Type Inhibitor

Formula $C_{18}H_{19}N_3O_3S$

MI 13: 8346

MW 357.43

Purity $\geq 99\%$

Solubility Soluble in ethanol and in buffered aqueous solutions (pH of 2.3).



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