

Losartan . potassium salt

AT1 antagonist.

Ordering Information

[Order Online »](#)

LKT-L5873-G001	1g
LKT-L5873-G005	5g
LKT-L5873-G025	25g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

CAS 124750-99-8

Formula $C_{22}H_{22}ClN_6O \cdot K$

MW 461.01

Purity $\geq 98\%$

Solubility Soluble in water to 100 mM, DMSO to 100 mM, and alcohol.



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-
be@enzolifesciences.com](mailto:info-be@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)