

# ADU-S100 sodium salt

Synthetic cyclic dinucleotide agonist of STING.

## Ordering Information

[Order Online »](#)

LKT-A145823-M001	1mg
LKT-A145823-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

<b>Alternative Name</b>	ML RR-S2 CDA Sodium Salt, STING-Inducer-1, MIW815, GTPL10192, Dithio-(Rp,Rp)-[cyclic[A(2',5')pA(3',5')p]]
<b>CAS</b>	1638750-95-4
<b>Formula</b>	$C_{20}H_{22}N_{10}O_{10}P_2S_2 \cdot 2 Na$
<b>MW</b>	734.5
<b>Purity</b>	≥99%
<b>Solubility</b>	Soluble in water (>73 mg/mL, >100 mM)



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)