

# DT MicroLoops LD – Small Aperture Assortment

Not sold in the U.S. or Japan.

5 each of 20, 35, 50 and 75  $\mu\text{m}$  aperture loops, SPINE length

## Ordering Information

[Order Online »](#)

JBS-M5-L18SP-A1LD	20ea
-------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details



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)