

# Cryo-EM Puck Gen 2.0

## (1 pc.)

The Cryo-EM Pucks Gen 2.0 have enhanced features over other cryo-EM pucks on the market. These include: 12 numbered well locations with easy-access tweezer notches Magnets in the bottom of the pucks, which securely hold the puck to the shelves of the shipping and storage canes Front grabbing tong location 2D barcoding for advanced sample tracking

### Ordering Information

[Order Online »](#)

JBS-X-CEMG2-1	1ea
---------------	-----

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

### Technical Info / Product Notes

#### **Puck Color:**

Selectable from blue, black, brown, gold, green, grey, orange, purple, red, violet

**Please indicate the desired color(s) upon ordering.**

Standard for Cat. No. X-CEMG2-1 is blue

#### **Puck Engraving:**

Standard is 6 characters, e.g. ABC-123

**Please indicate the desired engraving upon ordering.**



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)