

# Uni-Puck

The Uni-Puck is a sample pin storage & shipping container compatible with many automated sample mounting systems. It holds 16 sample bases/pins and consists of two parts: the sample enclosure and the sample base piece. Each part of the Uni-Puck has a unique serial number for identification.

## **Puck Color:**

Selectable from blue, black, brown, gold, green, grey, orange, purple, red, magenta, rainbow (for a set of 7 Uni-Pucks): 1x blue, 1x black, 1x green, 1x grey, 1x orange, 1x purple, 1x red.

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

Standard color is black.

## **Puck Engraving:**

Standard is 6 characters, e.g. ABC-001

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

## **Puck 2D-Barcode:**

Optional, ***please indicate (y/n) with your order.***

## **Ordering Information**

[Order Online »](#)

JBS-X-CC-110	1ea
--------------	-----

## **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)