

# 5-Vinyl-2'- deoxyuridine (5-VdU) (10 mg)

## Ordering Information

[Order Online »](#)

JBS-CLK-050-10	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5-Ethenyl-2'-deoxyuridine, 2'-Deoxy-5-vinyluridine, 2'-Deoxy-5-ethenyluridine

**Appearance** white to off-white solid

**Application Notes** DNA synthesis monitoring

**CAS** 55520-67-7

**Formula**  $C_{11}H_{14}N_2O_5$

**MW** 254.24

**Purity** ≥ 98 % (HPLC)

**Solubility** DMSO, methanol

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)