

5-Vinyl-2'- deoxyuridine (5-VdU) (50 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-050-50	50mg
----------------	------

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com