

5-FAM-Alkyne (5 x 5 mg)

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| JBS-CLK-057-25 | 5x5mg |
|----------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|---------------------|
| Handling | Store dark. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|--------------------|---|
| Appearance | yellow solid |
| CAS | 510758-19-7 |
| Emission Maximum | 513 nm |
| Excitation Maximum | 490 nm |
| Formula | $C_{24}H_{15}NO_6$ |
| MW | 413.38 |
| Purity | > 95 % (HPLC-MS, H NMR) |
| Solubility | aqueous buffer (pH >8), alcohols, DMSO, DMF |

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