

# Despropionylfentanyl

A major urinary metabolite of fentanyl

Major urinary metabolite of the narcotic analgesic fentanyl.

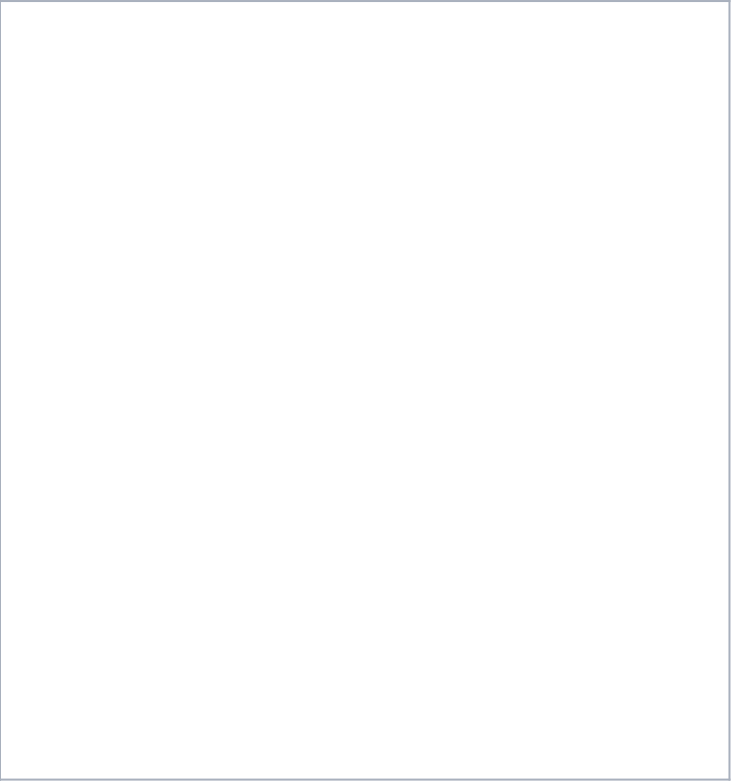
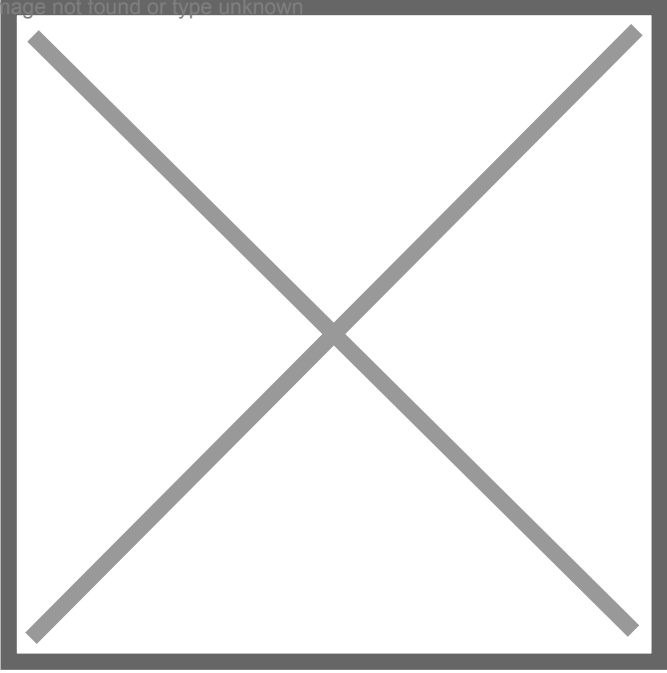
Citations: 1 [View Online »](#)

Ordering Information [Order Online »](#)

|               |      |
|---------------|------|
| BML-J110-0001 | 1mg  |
| BML-J110-0010 | 10mg |

Manuals, SDS & CofA [View Online »](#)

Image not found or type unknown



## Handling & Storage

|                   |   |
|-------------------|---|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months. |
| Long Term Storage | Ambient   |
| Shipping          | Ambient Temperature   |

**Regulatory Status** RUO - Research Use Only

## Product Details

|            |                          |
|------------|--------------------------|
| Appearance | Off-white solid.         |
| CAS        | 21409-26-7               |
| Formula    | $C_{19}H_{24}N_2$        |
| MW         | 280.4                    |
| Purity     | ≥99% (HPLC)              |
| Solubility | Soluble in 100% ethanol. |
| Source     | Synthetic.               |

Last modified: May 29, 2024

Image not found or type not known  
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)