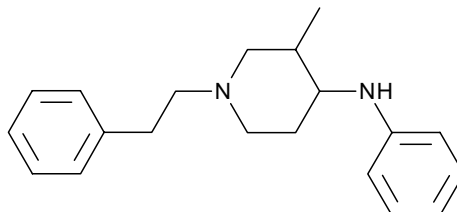


PRODUCT: Despropionyl-3-methylfentanyl

CATALOG NO.: J-115

LOT#: temp

STRUCTURE:



PHYSICAL APPEARANCE: Tan waxy solid (mp = 48-53°C)

MOLECULAR FORMULA: C₂₀H₂₆N₂

MOLECULAR WEIGHT: 294.4

PURITY: 99% (TLC) ; 50/50 mixture of *cis/trans* diastereomers

SOLUBILITY: Soluble in ethanol

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Major urinary metabolite of 3-methylfentanyl.¹

REFERENCES: 1. W.R.Hammargren and G.L.Henderson *J.Anal.Toxicol.* 1988
12 183

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.