

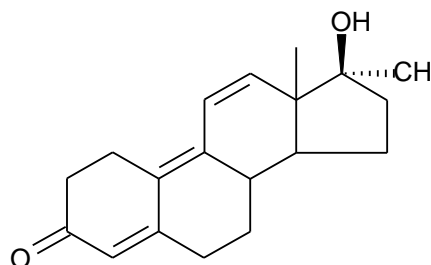
Methyltrienolone (R1881)

Product Number: NLP005005MG

Packaging: Amber screw-cap vial.

Storage Recommendations:

Methyltrienolone stored at -20°C away from light



MW. 284.38

Chemical Purity:

1. Thin-layer chromatography on silica gel using:

toluene : ethyl acetate, (8:2)

To stock purity is determined to be >97%, by UV visualization of one peak at 254nm.

2. NMR analysis performed for compound identification.

Special Instructions:

To increase the solubility of this product in aqueous solutions, pre-dissolve in a minimal amount of ethanol or DMSO.

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

NLP005005MG-R REV01