

# PRODUCT DATA SHEET



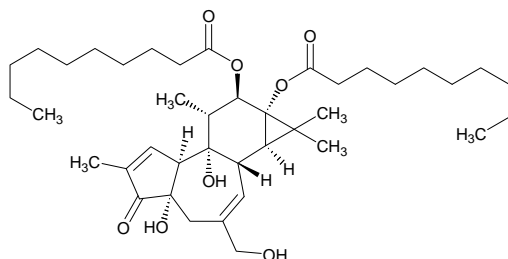
PRODUCT: 4 $\alpha$ -Phorbol 12,13-didecanoate

CAS NO: 27536-56-7

CATALOG NO.: PE-147

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Colorless oil

MOLECULAR FORMULA: C<sub>40</sub>H<sub>64</sub>O<sub>8</sub>

MOLECULAR WEIGHT: 672.9

PURITY: 99%

SOLUBILITY: Soluble in DMSO

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

APPLICATION NOTES: Use caution – product is a vesicant. Negative control for phorbol 12,13-didecanoate.<sup>1</sup>

REFERENCES: 1. M.Castagna *et al. J.Biol.Chem.* 1982 **257** 7847

**CAUTION!!!** Avoid all forms of contact with this product by wearing appropriate protective clothing (gloves, mask, lab coat). 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. The Material Safety Data Sheet (MSDS), is available on request.