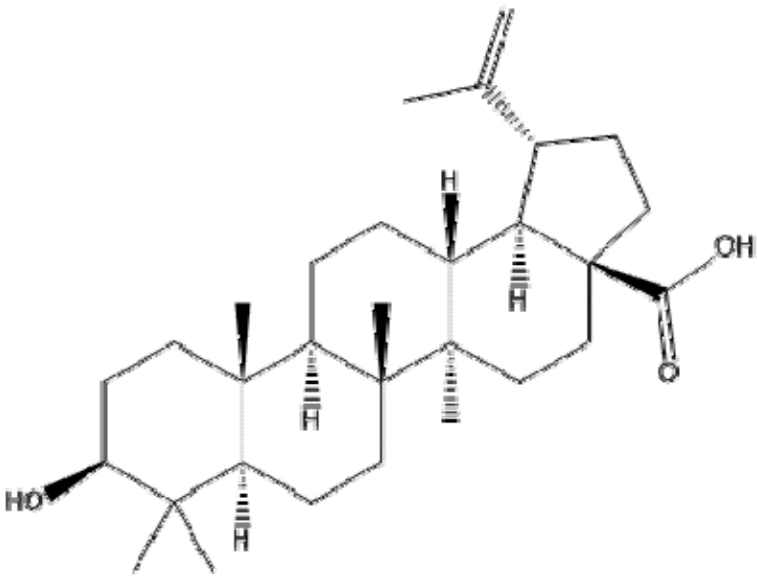


## Instruction Manual

Catalog Number	PK-CA577-1552-25
Description	Anti-tumor and anti-HIV agent. Induces apoptosis. Also displays TGR5 agonist activity (EC <sub>50</sub> = 1.04 μM).
Quantity	25 mg
Sequence / Molecular Weight / Molecular Formula	456.7 Da; C <sub>30</sub> H <sub>48</sub> O <sub>3</sub>
	
Purity	≥ 95% as determined by HPLC.
Appearance / Formulation / Solubility	White to off-white powder. Soluble in DMSO.
Storage & Stability	Store powder desiccated at -20°C for up to one year.
Applications	see Description
References	Genet, C., <i>et al.</i> (2010). <i>J. Med. Chem.</i> <b>53</b> , 178-190.
Caution	NA

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.