

(2Z)-4-[(1R,3aS,5S,12aS)-9-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-3a,4,5,7-tetrahydro-8,10-dihydroxy-3,3-dimethyl-11-(3-methyl-2-buten-1-yl)-7,13-dioxo-1,5-methano-1H,3H-furo[3,4-d]xanthen-1-yl]-2-methyl-2-butenoic acid

Expiration date: 29NOV2029

Results

Conforms

Reviewed and approved by: Ryan Seiffert



WWW.CAYMANCHEM.COM