

5Z,8Z,11Z,14Z-eicosatetraenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1Z)-1-octadecen-1-yloxy]ethyl ester

Expiration date: 24FEB2028

Tests

Results

Conforms

Reviewed and approved by: Ryan Seiffert



Copyright Cayman Chemical Company, 10/12/2018

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM