

2,2'-((2-(2-(2-(bis(carboxylatomethyl)amino)-5-(3-carboxylato-4-(2,7-dichloro-6-oxido-3-oxo-3H-xanthen-9-yl)benzamido)phenoxy)ethoxy)phenyl)azanediyldiacetate, hexapotassium salt

Purity Specification:	≥90%				
Molecular formula:	C43H27Cl2N3O16 • 6K				
CAS number:	Formula weight:	1147.20	Expiration date:	03APR2029	

Tests	Results
HPLC	Purity: 96.0 %

Reviewed and approved by: Michelle Riley

MSK

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2018

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM