



Calcium Detection Reagents

Product	Size	Catalog #
Calcein	100 mg	89200
Calcein, AM	1 mg	89201
Calcein, AM *UltraPure Grade*	1 mg	89202
Calcein, AM *UltraPure Grade* *Small Package*	50 µg x 20	89204
Calcein, AM *UltraPure Grade* *5 mM solution in anhydrous DMSO*	200 µL	89203
Calcein blue, AM	1 mg	89205
Coelenterazine *UltraPure Grade*	0.25 mg	82255
Coelenterazine cp *UltraPure Grade*	0.25 mg	82256
Coelenterazine f *UltraPure Grade*	0.25 mg	82257
Coelenterazine h *UltraPure Grade*	0.25 mg	82258
Coelenterazine hcp *UltraPure Grade*	0.25 mg	82259
Coelenterazine n *UltraPure Grade*	0.25 mg	82260
Dihydrocalcein, AM	1 mg	85700
Dihydrocalcein, AM *Small Package*	50 µg x 20	85701
Fluo-3, AM	1 mg	84024
Fluo-3, AM *UltraPure Grade*	1 mg	84025
Fluo-3, AM *UltraPure Grade* *Small Package*	50 µg x 20	84026
Fluo-3, AM *UltraPure Grade* *1 mM solution in anhydrous DMSO*	1 mL	84027
Fluo-3, pentaammonium salt	1 mg	84020
Fluo-3, pentapotassium salt	1 mg	84021
Fluo-3, pentasodium salt	1 mg	84022
Fura-2, AM	1 mg	84015
Fura-2, AM *UltraPure Grade*	1 mg	84016
Fura-2, AM *UltraPure Grade* *Small Package*	50 µg x 20	84017
Fura-2, AM *UltraPure Grade* *1 mM solution in anhydrous DMSO*	1 mL	84018
Fura-2, pentapotassium salt	1 mg	84011
Fura-2, pentasodium salt	1 mg	84012
Indo-1, AM	1 mg	84006
Indo-1, AM *UltraPure Grade*	1 mg	84007
Indo-1, AM *UltraPure Grade* *Small Package*	50 µg x 20	84008
Indo-1, AM *UltraPure Grade* *1 mM solution in anhydrous DMSO*	1 mg	84009
Indo-1, pentapotassium salt	1 mg	84000
Indo-1, pentasodium salt	1 mg	84001
Rhod-2, AM	1 mg	84034
Rhod-2, AM *UltraPure Grade*	1 mg	84035
Rhod-2, AM *UltraPure Grade* *Small Package*	50 µg x 20	84036
Rhod-2, AM *UltraPure Grade* *1 mM solution in anhydrous DMSO*	1 mL	84037
Rhod-2, tripotassium salt	1 mg	84030
Rhod-2, trisodium salt	1 mg	84031

Calcium detection reagents are also available in bulk packaging; please contact service@anaspec.com for more information.