

## FLIPR Tetra Instrument LED and Filter Combinations

Target Assay	LED (nm)	LED Part #	Emission Filter (nm)	Tetra Filter Part #	Tetra Kit Part #	Old Tetra Filter Part #	Old Tetra Kit Part #
Fura-2	335-345 and 380-390 (UV)	0200-6272 0200-6273	475-535	0200-6211	0200-6271	0200-6149	0200-6270
MQAE	360-380 (UV)	0200-6178	400-460	0200-6213		0200-6179	
Voltage Sensor Probe (VSP)	390-420 (UV)	0200-6135	440-480	0200-6205	0200-6208	0200-6136	0200-6137
Tango™ GPCR Assay System (FRET)	390-420 (UV)	0200-6135	440-480 515-575	0200-6205		0200-6136	
CFP, eCFP/YFP	420-455	0200-6148	475-535	0200-6211		0200-6149	
FLIPR Calcium Assay Kits	470-495 (Blue)	0200-6128	515-575	0200-6203	0200-6206	0200-6108	0200-6130
Calcium, Fluo-4	470-495 (Blue)	0200-6128	515-575	0200-6203	0200-6206	0200-6108	0200-6130
Calcium, Fluo-3	495-505	0200-6175	526-586	0200-6212		0200-6176	
JC-1	495-505	0200-6175	565-625	0200-6212		0200-6109	
FLIPR Membrane Potential Assay Kits	510-545 (Green)	0200-6127	565-625	0200-6204	0200-6207	0200-6109	0200-6129
Rhod-2 & Rhod-4	510-545	0200-6127	565-625	0200-6204		0200-6109	
Alexa 633 & Bodipy	610-626	0200-6150	646-706	0200-6214		0200-6170	
BRET 2 luminescence assay with ICCD Camera	NA	NA	440-480 526-586	0200-6205 0200-6212		NA NA	
Single Custom Filter Holder	NA	NA		0200-6276		NA	
Three Custom filter holders	NA	NA		0200-6221		0200-6119	