



FLIPR® Penta and Tetra Systems

Instrument LED and filter combinations

Target Assay	LED (nm)	Penta and Tetra LED Part #	Emission Filter (nm)	Penta Filter Part #	Penta Kit Part #	*Tetra Filter Part #	*Tetra Kit Part #
Fura-2 QBT Calcium Kit	335-345 and 380-390 (UV)	5201196 0200-6273	475-535	0200-6211	5302739	0200-6149	0200-6270
MQAE	360-380 (UV)	0200-6178				0200-6179	
Voltage Sensor Probe (VSP)	390-420 (UV)	0200-6135	440-480 565-625	0200-6205 0200-6204	0200-6208	0200-6136 0200-6109	0200-6137
Tango™ GPCR Assay System (FRET)	390-420 (UV)	0200-6135	440-480 515-575	0200-6205 0200-6203		0200-6136 0200-6108	
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	0200-6204		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 Luminescence Assay with HSEMCCD Camera	NA	NA	440-480 526-586	0200-6205 0200-6212		NA NA	
Three custom filter holders	NA	NA		0200-6221		0200-6119	

*Tetra filter part and kit part for instruments with serial number less than 200. For Tetra FT0200 + (include 200), use the Penta EM filter part numbers.

Emission filters specifications for 0200-6221 FLIPR Custom Filter Kit

Suggested vendors

Omega Filters

Kim Cook

Delta Campus, 21 Omega Drive, Brattleboro, VT 05310

+1.866.488.1064 x8313 (work) | +1.802.251.7313 (mobile) | kcook@omegafilters.com

Semrock

Sabine Blessing, OEM Account Manager

3625 Buffalo Road, Suite 6, Rochester, NY 14624

+49.1590.4318499 (mobile) | sblessing@idexcorp.com

Specifications

- UV to 1.2 micron blocking: average OD = 5 (block stray light)
- Optical polishing and grinding
- Less than 40 arcsec parallelism
- Anti-reflection coating (AR) both sides
- 2.441" +.000/-.010 diameter (62.0 mm, clear aperture 58 mm)
- Maximum filter thickness .275" (6.98 mm)

Additional notes

- Recommended minimum average transmission in the passband: 75%
- There is a possibility of light around 800–900 nm that needs to be blocked, so avoid using a passband that goes up to these wavelengths.
- Typical lead time is two weeks for filters which are already made (previous run); or six weeks for custom filters.



Contact Us

Phone: +1.800.635.5577

Web: www.moleculardevices.com

Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1.800.635.5577

United Kingdom +44.118.944.8000

Europe* 00800.665.32860

China +86.4008203586

Taiwan/Hong Kong +886.2.2656.7585

Japan +81.3.6362.9109

South Korea +82.2.3471.9531

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland