

We've got your wavelength covered



Mix and match to detect your favorite assays on the FLIPR® System

FLIPR Tetra Instrument configurations

	Target assay application	Ex LED (nm)	Em filter (nm)
Fluorescence	New! Fura-2	335-345 and 380-390	475-535
	MQAE	360-380	400-460
	Voltage Sensor Probe (VSP)	390-420	440-480
	Tango GPCR assay system (FRET)	390-420	440-480, 515-575
	CFP; eCFP/YFP	420-455	475-535
	FLIPR® Calcium Assay	470-495	515-575
	Calcium (Fluo-4), Calcium (Fluo-8)	470-495	515-575
	Calcium (Fluo-3)	495-505	526-586
	JC-1	495-505	565-625
	FLIPR® Membrane Potential Assay	510-545	565-625
	Rhodamine-2, Rhodamine-4	510-545	565-625
	Alexa 633 and Bodipy	610-626	646-706
Luminescence	Aequorin	Luminescence	
	GloSensor	Luminescence	
	BRET ²	Luminescence	440-480, 526-586

We offer a broad range of optics for the FLIPR® Tetra System to enable the application flexibility you require to support your research and screening. LEDs and filters are easily changed on the system in just a few minutes. Move across the ultraviolet and visible light spectra, to luminescence, and back with no requirement for a service visit.

Scan the QR code to request more information or visit www.moleculardevices.com.



Additional dyes and applications may be addressed. Ask us for more information about your applications of interest. Custom filter holders are available to work with additional emission filters for your applications.

