

BioLegend Flow Cytometry Fluorophore Guide

Fluorophore	Emission Max	Similar Fluorophores	Brightness Index
Spark UV™ 387	387 nm	BUV395™	●●○○○
Brilliant Violet 421™	421 nm	Pacific Blue™, V450, eFluor® 450, AF405	●●●●●
Spark Violet™ 423	415 nm, 423 nm	Super Bright 436	●●●○○
Pacific Blue™	455 nm	BV421™, V450, eFluor® 450, AF405	●○○○○
Brilliant Violet 510™	510 nm	V500, AmCyan, Pacific Orange™	●●○○○
Spark Violet™ 538	538 nm	Pacific Orange™, Krome Orange™	●○○○○
Brilliant Violet 570™	570 nm	-	●●○○○
Brilliant Violet 605™	603 nm	Qdot 605	●●●●○
Brilliant Violet 650™	645 nm	Qdot 655	●●●●○
Brilliant Violet 711™	711 nm	Qdot 705	●●●●●
Brilliant Violet 750™	750 nm	-	●●●○○
Brilliant Violet 785™	785 nm	Qdot 800	●●●●○
Alexa Fluor® 488	519 nm	FITC, BD Horizon Brilliant™ Blue 515	●●●○○
KIRAVIA Blue 520™	520 nm	FITC, AF488, BD Horizon Brilliant™ Blue 515	●●●○○
FITC	525 nm	AF488, BD Horizon Brilliant™ Blue 515	●●●○○
Spark Blue™ 550	540 nm	AF532	●○○○○
Spark Blue™ 574	574 nm	-	●○○○○
PerCP	675 nm	-	●○○○○
PerCP/Cyanine5.5	690 nm	PerCP-eFluor™ 710, BD Horizon Brilliant™ Blue 700	●●○○○
PE	575 nm	-	●●●●●
PE/Dazzle™ 594	610 nm	BD Horizon™ PE-CF594, ECD, PE-Texas Red®, AF594	●●●●●
PE/Fire™ 640	639 nm	-	●●●●○
PE/Cyanine5	670 nm	-	●●●●●
PE/Fire™ 700	695 nm	PE/Cyanine5.5	●●●●●
PE/Cyanine7	774 nm	-	●●●●○
PE/Fire™ 810	806 nm	-	●●●●○
Spark YG™ 581	581 nm	cFluor® YG584	●●○○○
Spark YG™ 593	593 nm	BD Horizon RealYellow™ 586	●●○○○
APC	660 nm	AF647, eFluor™ 660	●●●●○
Alexa Fluor® 647	668 nm	APC, eFluor™ 660	●●●●○
Spark NIR™ 685	685 nm	AF660	●●●○○
Spark Red™ 718	711 nm	AF700, BD Horizon™ Red 718, APC-R700	●●●●○
Alexa Fluor® 700	719 nm	Spark Red™ 718, BD Horizon™ Red 718, APC-R700	●●○○○
APC/Cyanine7	774 nm	APC-H7, APC-eFluor™ 780	●●○○○
APC/Fire™ 750	774 nm	APC/Cyanine7, APC-H7, APC-eFluor™ 780	●●○○○
APC/Fire 810™	807 nm	-	●●○○○

Excitation Lasers				
Ultraviolet	Violet	Blue	Yellow/Green	Red

Viability Dyes	Emission Max	Similar Viability Dyes
Zombie UV™	459 nm	LIVE/DEAD™ Fixable Blue
Zombie Violet™	423 nm	LIVE/DEAD™ Fixable Violet
Helix NP™ Blue	470 nm	SYTOX™ Blue
Zombie Aqua™	516 nm	LIVE/DEAD™ Fixable Aqua
Zombie Yellow™	572 nm	LIVE/DEAD™ Fixable Yellow
Zombie Green™	515 nm	LIVE/DEAD™ Fixable Green
Helix NP™ Green	519 nm	SYTOX™ Green
Apotracker™ Green	520 nm	-
Zombie Red™	624 nm	LIVE/DEAD™ Fixable Red
Helix NP™ NIR	660 nm	SYTOX™ Red
Zombie NIR™	746 nm	LIVE/DEAD™ Fixable Far Red

BioLegend is the leading provider of flow cytometry reagents, bolstered by an extensive antibody clone library and spectrum-spanning fluorophores. Utilize our charts to build the perfect multicolor panel.

Learn more and see our updated fluor list at:
biolegend.com/en-us/fluorophore-families



05-0116-00



BioLegend is a registered trademark of BioLegend, Inc. All other trademarks are the property of their respective owners. AF=Alexa Fluor®, BV=Brilliant Violet™, BUV=Brilliant Ultraviolet™. For trademarks, visit: biolegend.com/en-us/trademarks