

VersaFluor™ Filter Selection Guide

Application	Excitation Filters	Emission Filters
DNA Quant (Hoechst 33258)	EX 360/40 (340–380 nm)	EM 460/10 (455–465 nm)
β-Galactosidase (MUG)	EX 360/40 (340–380 nm)	EM 460/10 (455–465 nm)
β-Glucuronidase (MUGluc)	EX 360/40 (340–380 nm)	EM 460/10 (455–465 nm)
Apopain (Z-DEVD-AFC)	EX 360/40 (340–380 nm)	EM 520/10 (515–525 nm)
pH indicator (BCECF)	EX 490/10 (485–495 nm)	EM 520/10 (515–525 nm)
Calcium indicator (Fluo-3)	EX 490/10 (485–495 nm)	EM 520/10 (515–525 nm)
DNA Quant (PicoGreen™)	EX 480/20 (470–490 nm)	EM 520/10 (515–525 nm)
Oligo Quant (OliGreen™)	EX 480/20 (470–490 nm)	EM 520/10 (515–525 nm)
Protein Quant (NanoOrange™)	EX 480/20 (470–490 nm)	EM 590/10 (585–595 nm)
Fluorescein	EX 490/10 (485–495 nm)	EM 520/10 (515–525 nm)
Ethidium Bromide	EX 510/10 (505–515 nm)	EM 620/10 (615–625 nm)
RNA Quant (RiboGreen™)	EX 490/10 (485–495 nm)	EM 520/10 (515–525 nm)