



Bigfoot Cell Sorter instrument configuration

Choose your configuration of up to 9 lasers and 60 channels

Instrument configurations offering spectral analysis and conventional compensation range from 5 to 9 lasers and up to 60 parameters, while configurations offering conventional compensation include 4 or 5 laser models and up to 35 parameters. Apart from lasers and detectors, all instrument configurations share the same technical specifications.



Invitrogen™ Bigfoot™ Cell Sorter configurations

All spectrally enabled instrument configurations also offer conventional compensation capabilities.

Number of fluorescence detectors	4-laser conventional (Cat. No. PL00304)	5-laser conventional (Cat. No. PL00303)	5-laser spectral (Cat. No. PL00302)	6-laser spectral 785 nm (Cat. No. PL00301)	6-laser spectral 445 nm (Cat. No. PL00300)	7-laser spectral (Cat. No. PL00299)	9-laser spectral (Cat. No. PL00285)
348 nm	387/11		•	•	•	•	•
	420/10		•	•	•	•	•
	434/17			•	•	•	•
	455/14			•	•	•	•
	465/22		•		•	•	•
	473/15			•	•	•	•
	507/19			•	•	•	•
	525/35		•				
	549/15			•	•	•	•
	575/15			•	•	•	•
	583/30		•				
	615/24			•	•	•	•
	670/30		•	•	•	•	•
	690LP		•				
	728/40			•	•	•	•
	750LP			•	•	•	•
405 nm	420/10	•	•	•	•	•	•
	434/17			•	•	•	•
	455/14			•	•	•	•
	465/22	•	•				
	473/15			•	•	•	•
	507/19			•	•	•	•
	525/35	•	•				
	549/15			•	•	•	•
	575/15			•	•	•	•
	583/30	•	•				
	615/24			•	•	•	•
	661/20		•				
	710/20			•	•	•	•
	720/60	•	•				
	747/33			•	•	•	•
	750LP	•	•				
770LP			•	•	•	•	
445 nm	465/22				•	•	•
	525/35				•	•	•
	583/30					•	•
	650LP					•	•
488 nm/632 nm (BL)	507/19	•	•	•	•	•	•
	549/15	•	•	•	•	•	•
	583/30	•	•	•	•	•	•
	593/52		•				
	615/24	•		•	•	•	•
	670/30	•		•	•	•	•
	692/80		•				
	720/60	•		•	•	•	•
561 nm/594 nm (BL)	750LP	•	•	•	•	•	•
	575/15	•	•	•	•	•	•
	589/15	•	•	•	•	•	•
	605/15			•	•	•	•
	615/24	•	•				
	622/15			•	•	•	•
	661/20	•	•				
	685/15			•	•	•	•
	700/13	•	•				
	720/24	•	•				
	750LP	•	•				
	760/50			•	•	•	•
	800/12			•	•	•	•
	832/37			•	•	•	•
640 nm	860LP			•	•	•	•
	670/30	•	•	•	•	•	•
	700/13	•	•	•	•	•	•
	720/24			•	•	•	•
	728/40		•				
	750LP	•	•				
785 nm	760/50			•	•	•	•
	770LP			•	•	•	•
	800/12			•	•	•	•
	832/37			•	•	•	•
860LP			•	•	•	•	

Learn more at thermofisher.com/bigfoot

