

Brilliant Violet™ 421

DAPI

*Pacific Blue™

*Hoechst

*Alexa Fluor® 488/FITC

Pacific Orange™

Cy™3

Alexa Fluor® 555

*PE

Texas Red®

PE-Texas Red®

*Alexa Fluor® 594

*PI/7-AAD

APC

Max. Emission Wavelength (nm)

421

Max. Excitation Wavelength (nm)

407

Excitation Laser Line (nm)

405

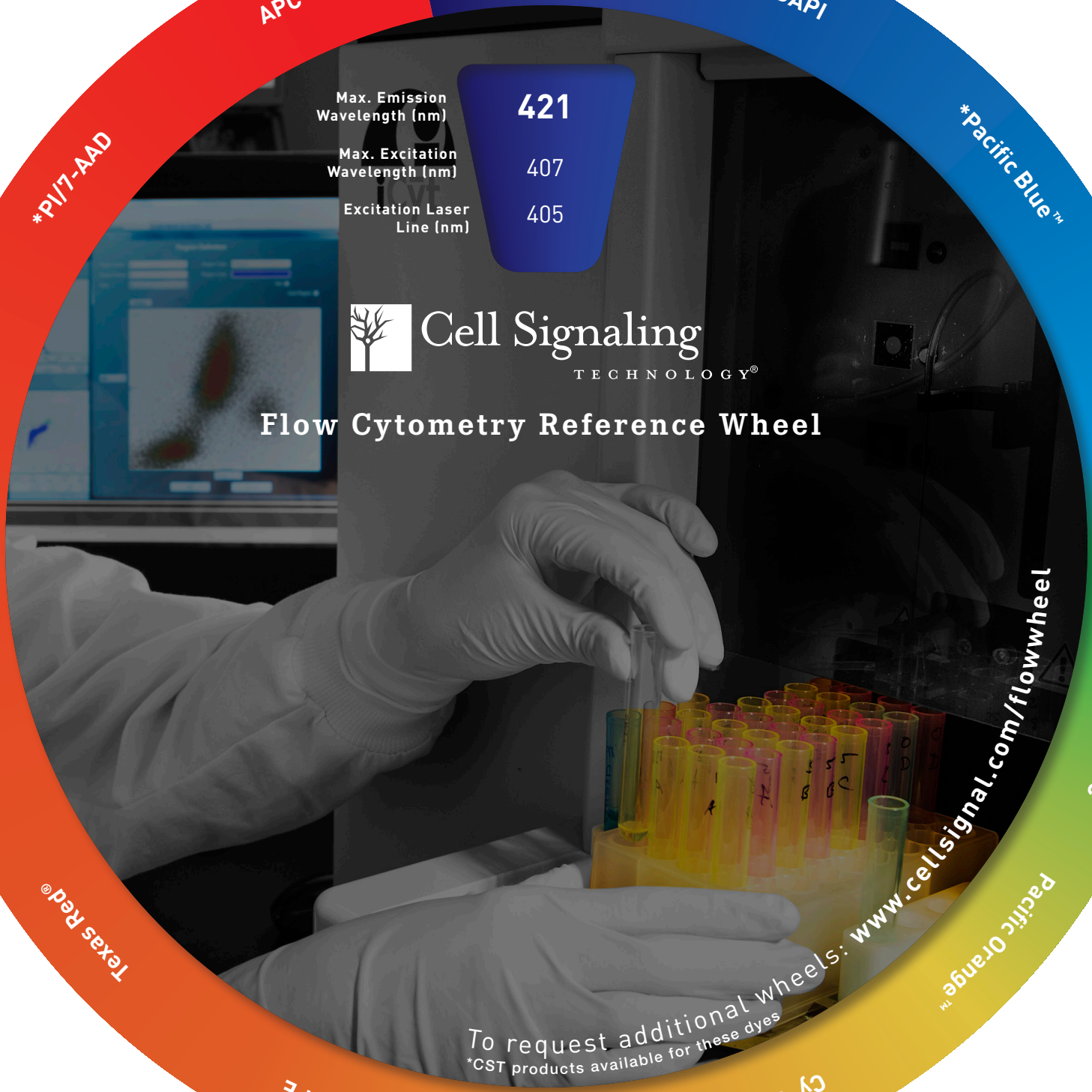


Cell Signaling

TECHNOLOGY®

Flow Cytometry Reference Wheel

To request additional wheels: www.cellsignal.com/flowwheel
*CST products available for these dyes



Alexa Fluor® 647*

PE-Cy™5

PerCP

*DRAQ7™

*DRAQ5®

PE-Cy™5.5

APC-Cy™5.5

PerCP-Cy™5.5

*DyLight™ 680

Alexa Fluor® 700

APC-Cy™7

PE-Cy™7

*DyLight™ 800

Alexa Fluor® 790

Max. Emission Wavelength (nm)

665

Max. Excitation Wavelength (nm)

650

Excitation Laser Line (nm)

633/635

Tips for selecting fluorochromes for multicolor flow cytometry

Research the fluorochrome properties – ex. tandem dyes are sensitive to photobleaching and can become uncoupled; some fluorochromes are better suited for intracellular staining than others (ex. PE-Cy™5 can withstand the harsh washing conditions required for intracellular staining).

To limit the amount of compensation required, select a single fluorochrome in each laser excitation range, rather than exciting multiple fluorochromes with the same laser.

Choose fluorochromes with the least amount of spectral overlap.

Select the brightest fluorochromes that can be detected by your instrument.

Pair the brightest fluorochromes with the lower expressing proteins and the dimmer fluorochromes with the highly expressed proteins.



Cell Signaling

TECHNOLOGY®

For additional flow cytometry resources: www.cellsignal.com/flow

5W000033P

Cell Signaling Technology® is a trademark of Cell Signaling Technology, Inc. / Alexa Fluor®, Texas Red®, Pacific Blue™, and Pacific Orange™ are trademarks of Molecular Probes. / Cy™ is a trademark of GE Healthcare. / Brilliant Violet™ is a trademark of Becton, Dickinson and Company. / DRAQ5® and DRAQ7™ are trademarks of Bioss Limited. / DyLight™ is a trademark of Thermo Fisher Scientific, Inc. and its subsidiaries.