

Discover what others can only imagine!

BD CellView™ Technology Applications



ANALYZE CELL-CELL INTERACTIONS IN REAL TIME



Gain Deeper Immunological Insights

BD CellView™ Technology, part of the BD FACSDiscover™ Platform, takes you beyond conventional flow cytometry analysis. By combining spectral flow cytometry with real-time imaging, visualize and sort immunological synapses as they form, unlocking deeper understanding, richer datasets and faster paths to publication.

Challenges you might be facing

Finding questionable population in flow cytometry like CD3+CD14+ double positive events?

- » **Are these artifacts or biology?**
- » **Does imaging enable the identification of these events?**

If you're studying immune interactions, every missed synapse is a missed insight, and potentially a missed discovery. Traditional flow cytometry often discards doublets as artifacts, making it nearly impossible to confidently identify and sort true cell-cell interactions. That means critical data about how T cells kill tumors, how macrophages polarize and how cancer evades immune attack could be slipping through your fingers.

Don't miss out on seeing what others can't

Here's what you gain with BD CellView™ Technology:

- » **See what others can't:** Precisely identify true doublets in real time.
- » **Reveal hidden complexity:** Visualize immune synapses using imaging parameter.
- » **Distinguish true biology from noise:** Combine label-free and fluorescence imaging to separate real synaptic events from coincident artifacts.

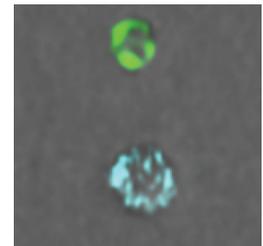
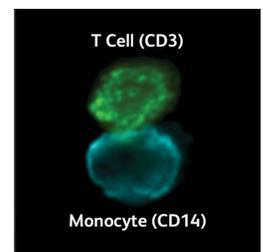
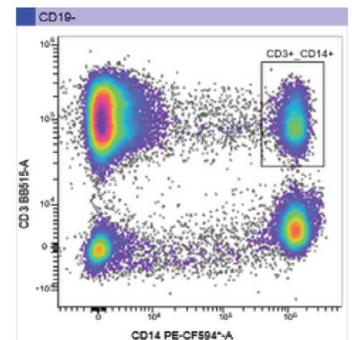
This technology in conjunction with BD FACSCorus™ Software, allows for real time assessment of your sample so you can visually confirm your gating decisions as you go. It allows you to confidently sort your population and ensure your data is accurately and truly representative of the biology you are trying to understand.

Don't only imagine! Discover how spatial context drives higher-dimensional analysis

- » **Detect true immunological synapses through integrated imaging and flow cytometry.**

Cut through the noise! By integrating spectral flow cytometry with real-time imaging, researchers can confidently distinguish true immunological synapses from coincident events, enabling precise characterization of cell-cell interactions critical for immuno-oncology.

Spatial meets spectral



- » **Reveal hidden events and see synapse interactions through image parameter information and make spotting overlapping synapses faster and more accurate.**

Automatically generated imaging parameters, such as correlation of fluorescent channels, uncover hidden or overlapping synapse interactions that traditional gating might miss, providing deeper insight into immune cell communication.

- » **Don't miss on critical synaptic events. Identify synaptic events without the need for labels or dyes.**

See more biology in its purest form. The BD FACSDiscover™ Instruments enable researchers to reveal critical synaptic events without relying on labels or dyes, using advanced image parameters for confident identification. This label-free approach allows scientists to observe immune cell interactions in their purest biological form, preserving native complexity while accelerating discovery.

From insight to impact: stay ahead of the curve with BD CellView™ Technology

If you rely on traditional flow cytometry, you might be missing critical insights hidden in complex cell-cell interactions, insights that could fuel your next publication or even your next big discovery. BD CellView™ Image Technology with BD FACSDiscover™ S8 Cell Sorter and BD FACSDiscover™ A8 Cell Analyzer, puts you at the forefront of innovation by integrating spectral flow cytometry with spatial and morphological detail.

Stay ahead, publish faster and accelerate your career with technology designed for the next generation of immunology research.

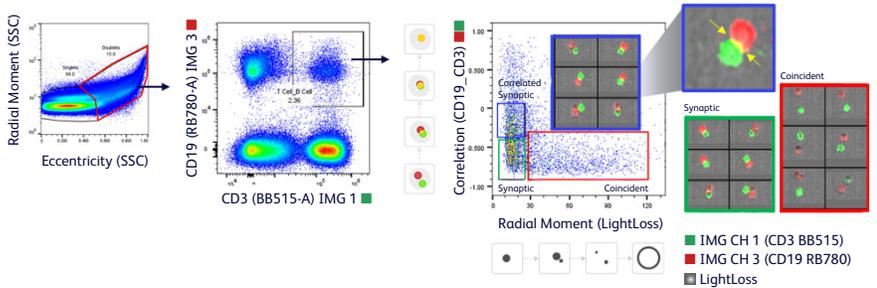
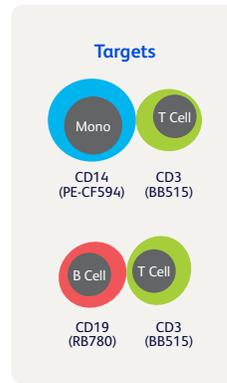
Check the following resources to learn more

<https://www.bdbiosciences.com/cellview>

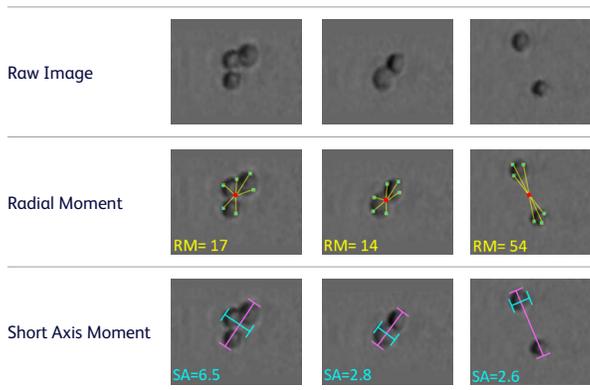
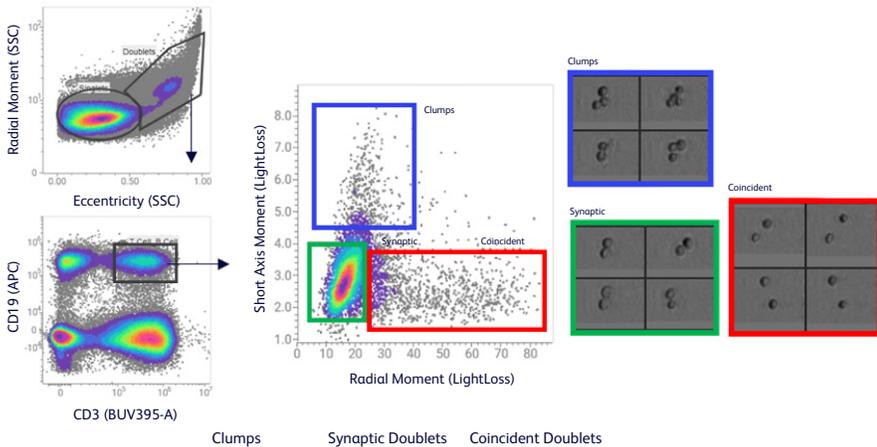
<https://www.bdbiosciences.com/en-us/learn/applications/cell-view-image-technology?tab=videos>

Cell-cell interaction: combining immunophenotyping and imaging

Cell Interaction: 20 Color Panel		
Laser	Antigen	Fluorochromes
UV	CD4	BUV395
	CD27	BUV563
	CD28	BUV615
	TCR $\gamma\delta$	BUV661
	CD39	BUV737
Violet	CD8	BUV805
	CD197	BV421
	CD45	BV480
	CD16	BV605
	IgG	BV750
Blue	CD45RA	BV786
	CD3	BB515
	CD14	PE-CF594
Yellow/Green	CD19	RB780
	CD25	RY586
Red	CD20	RY780
	CD127	AF647
	IgM	R718
	HLA-DR	APC-H7



Cell-cell interaction: Label free analysis



BD flow cytometers are Class 1 Laser Products. For Research Use Only. Not for use in diagnostic or therapeutic procedures.

BD Biosciences, Milpitas, CA 95035, USA | bdbiosciences.com

BD, the BD Logo, BD CellView, BD FACSDiscover and BD SpectralFX are trademarks of Becton, Dickinson and Company or its affiliates. © 2025 BD. All rights reserved. BD-167597 (v1.0) 0126

