

Discover valuable insights on breakthroughs, innovative strategies, and cell-based approaches for immunotherapeutic discovery and development.

Immunotherapeutics including immuno-oncology (IO) drugs is a class of treatment modality designed to augment the body's natural immune defense against cancer. This field is currently dominated by therapeutics development programs targeting immune checkpoint modulators and cytokines as these molecules hold significant promise in cancer therapy. Another key modality is the modulation of Fc-mediated effector mechanisms of therapeutic antibodies, including antibody-dependent cell-mediated cytotoxicity (ADCC), antibody-dependent cell phagocytosis (ADCP), complement dependent cytotoxicity (CDC), bi-specific antibody-mediated T-cell redirection, chimeric antigen receptor T-cell (CAR-T), adoptive T-cell therapies, and other cytotoxicity-focused applications.

Learn how you can rapidly screen and develop small molecules, biologics (peptides, antibodies, or bispecific antibodies), and cellular therapies for your immunotherapeutic discovery and development program.

New eBook Benefits:

- Review the latest on immunotherapeutic evolution, innate and adaptive immune systems, regulating T-cell responses, harnessing cytokines, and role of cell-based assays
- Gain expert immunotherapeutic drug discovery and development insights via exclusive Q&A's from industry leader Dr. Paige Mahaney Sr. VP and Corporate Head of Drug Discovery Research at Exelixis
- Connect to curated articles and application notes on functional, cell-based assays for checkpoints, cytokines/interleukins, growth factor receptors, therapeutics mediating cellular cytotoxicity, and more
- Access case studies and white paper on immune checkpoint receptors and KILR® cytotoxicity bioassays
- Unlock knowledge-based videos and webinars on immunotherapeutic cell-based assays
- Explore resources on immunotherapeutic product types, applications, targets, and programs

Learn More at [→ discoverx.com](https://discoverx.com)



DOWNLOAD
New Immunotherapeutics eBook at
discoverx.com/immunother-ebook



Immuno-oncology



Cytotoxicity



Checkpoint Modulators



Cytokine/Interleukin Receptors