

CERTIFICATION PROGRAM FOR CELL-BASED ASSAYS

Take Your Bioassay Services to the Next Level



Biopharmaceutical companies developing and manufacturing biologics and biosimilars are in need of CRO/CDMO's that can support their therapeutic programs with robust bioassays. At Eurofins DiscoverX, we have the largest portfolio of qualified bioassays available and custom bioassay development services. With over two decades of experience, biopharma clients seek our advice when moving their drug product downstream and want a CRO/CDMO that can successfully run our bioassays for their therapeutic.

This Certification Program from Eurofins DiscoverX is designed to qualify CRO/CDMO's to run our bioassays. Certified CRO/CDMO's are then marketed to biopharma clients focused on their downstream therapeutic efforts.

Want to be able to offer the best bioassays to your clients? Read on to learn more.

VALUE FOR CERTIFIED CRO/CDMO's

- Pre-release of new assay targets before new products go into the Eurofins DiscoverX website
- Availability of Field Application Specialist (FAS) for additional onsite trouble-shooting or training (fees may apply)
- Access to generate exclusive MCB/WCB for bioassays that reduces the need for bridging studies
- Reservation of pre-qualified lots with a minimum of 6 months shelf-life
- Ability to place bulk orders for QC lot release testing clients



Begin Your Certification

1  **Contact us at**
CRO_Certification@discoverx.com

4  **Attend a 3-day theoretical and wet-lab workshop at your site**

2  **Schedule a workshop performed by Eurofins DiscoverX FAS team**

5  **Obtain your Certification**

3  **Sign agreement letter prior to attending workshop**

CRO = Contract Research Organization

CDMO = Contract Development and Manufacturing Organization

CERTIFICATION PROGRAM FOR CELL-BASED ASSAYS

REQUIREMENTS FROM APPLYING CRO/CDMO

- A commitment to incorporate Eurofins DiscoverX bioassays and cell lines into their assay development portfolio.
- Expectation to share application data with Eurofins DiscoverX (without violating client confidentiality) that could be used to show the value of these products.
- To co-sponsor seminars and publications as appropriate.
- A minimum annual purchase of cell-based assay products.

Illustrated 3-Day Theoretical and Wet-Lab Workshop

DAY 1



AM: Instructions and wet-lab prep
(bioassay 1 cell thawing and plating)



PM: Classroom discussions
(MOA, bioassay platform and protocols, and results analysis)

DAY 2



AM: Wet-lab
(bioassay 1 ligand addition and
bioassay 2 cell thawing and plating)



PM: Wet-lab
(bioassay 1 plate reading and
data analysis)

DAY 3



AM: Wet-lab
(bioassay 2 plate reading and data
analysis)



PM: Classroom
(troubleshooting assays and
completion of certification)

COMMITMENT FROM EUROFINS DISCOVERX

- Once you are a certified CRO/CDMO, when a client comes to us asking for recommendations for CRO/CDMO's experienced in our bioassays, we will provide your company's name/contact to the client (non-exclusive).
- Your name/logo will appear on our website and in our marketing materials, stating your company's status as a certified CRO/CDMO.
- You will have access to co-sponsor seminars, white-papers, and other customer-education materials.

Want to Get Started on Your Certification for Cell-Based Assay? Contact us at CRO_Certification@discoverx.com.