

Advances in tools to compare potency, efficacy, safety & quality of follow-on biologics or biosimilars

Nadia Tagnaouti, Ph. D., Helen Roberts, Ph. D., Abhi Saharia, Ph. D., and Jane Lamerdin, Ph. D.
DiscoverX Corporation, Fremont, CA 94538-3142

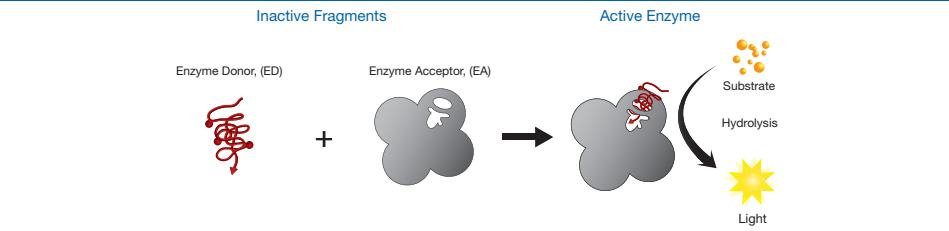
Abstract

Regulatory authorities define biosimilars as drugs that are “highly similar” to or “interchangeable” with an approved biologic. The lack of a pre-defined path to demonstrate similarity leaves the burden of proof on the biosimilar developers. Key requirements for efficacy, quality and potency testing often require a complex set of bioassays and/or cell-based assays that are also used to assess any unwanted clinical immune response to their biosimilar. This is particularly challenging to address for biosimilars.

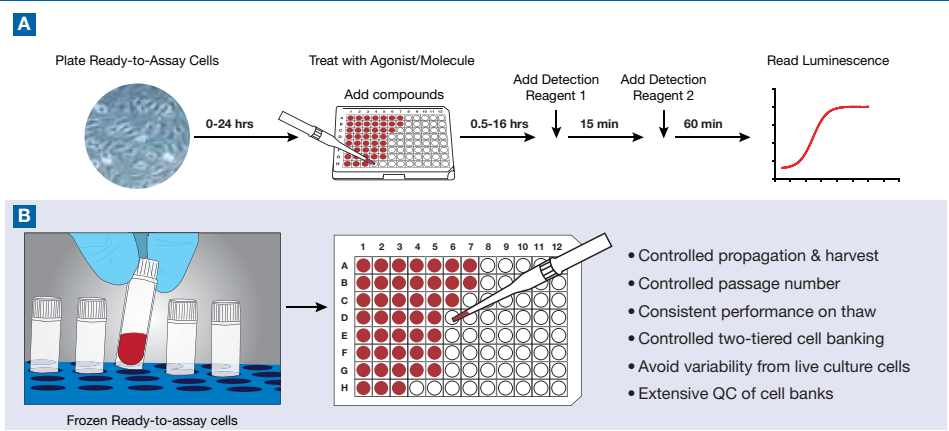
Commercially available cell-based assays for the development of biosimilars targeting different marketed drugs can significantly shorten development time while enabling easy adoption and greater reproducibility across multiple global sites. Examples discussed here include cell-based assays for Bevacizumab, highlighting how commercial tools can enable a biosimilar developer to move more quickly and be the first to market or submission.

Additionally, the emergence of complex human primary cell culture and co-culture *in vitro* models as well as biomarker-based phenotypic assays offer high-throughput and robust solutions for assessing mechanism of action and human safety and efficacy.

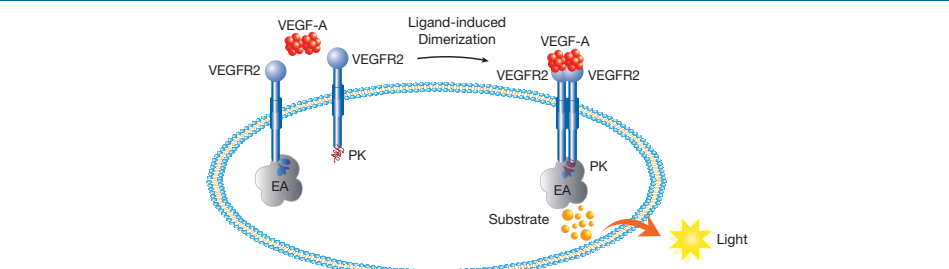
PathHunter Enzyme Fragment Complementation



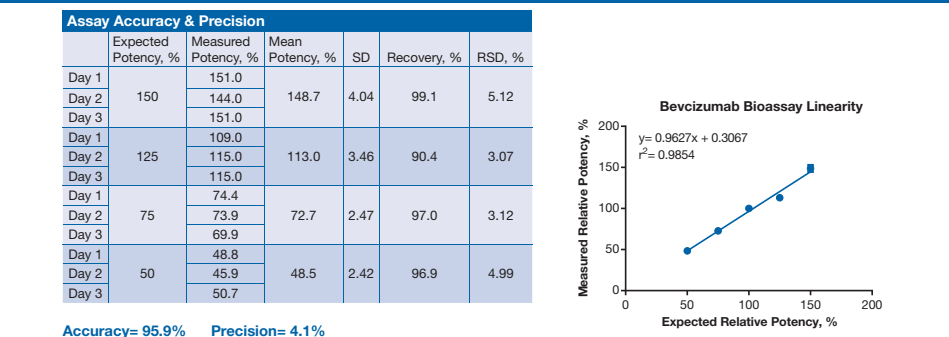
A Simple Homogenous Protocol With Rapid Results



PathHunter Bevacizumab Bioassay



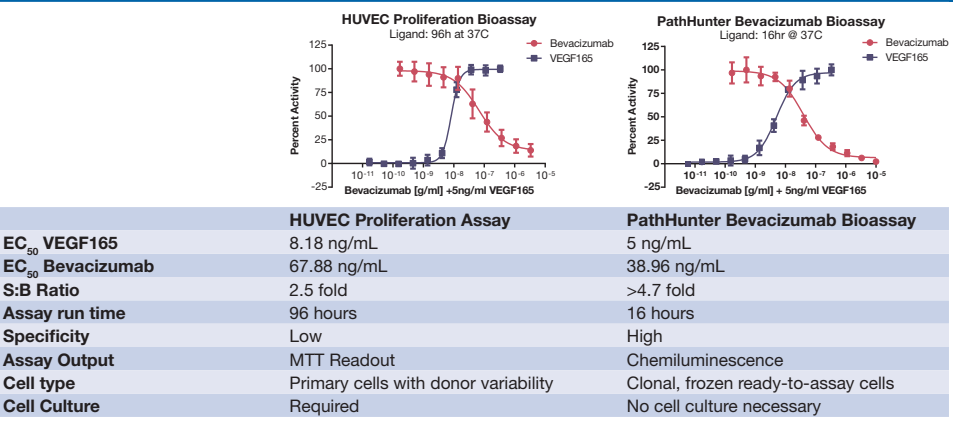
Excellent Accuracy and Precision for PathHunter Bevacizumab Bioassay



VEGF-A is known to cause homodimerization of VEGFR2 (KDR), as the first step in the activation cascade of these receptors. Anti-VEGF-A antibodies such as Avastin (Bevacizumab), Ranibizumab and Aflibercept bind to VEGF-A and prevent this dimer formation, leading to inhibition of VEGF-A dependent signaling. Here we have tested the VEGFR2 homodimer assay with VEGF-A, demonstrating a robust response and a high level of reproducibility with multiple runs.

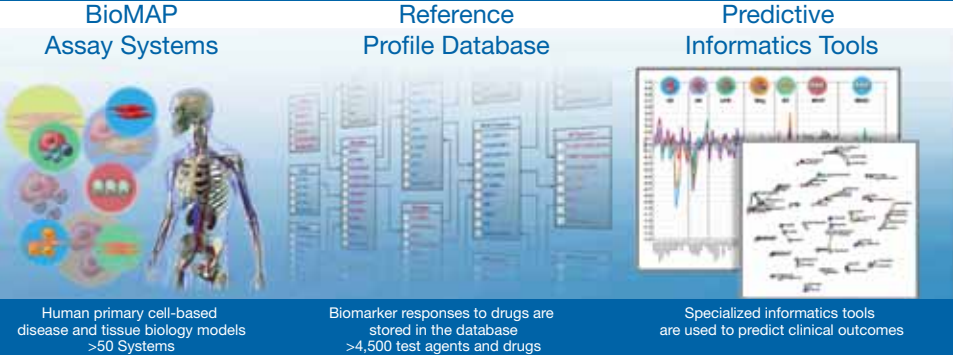
The VEGFR dimerization assay was tested with four test samples, from 50% to 150%, compared to a reference standard (100%). The measured relative potencies were plotted against the expected relative potencies with a very high degree of accuracy and precision.

HUVEC Proliferation Assay vs. PathHunter Bevacizumab Bioassay



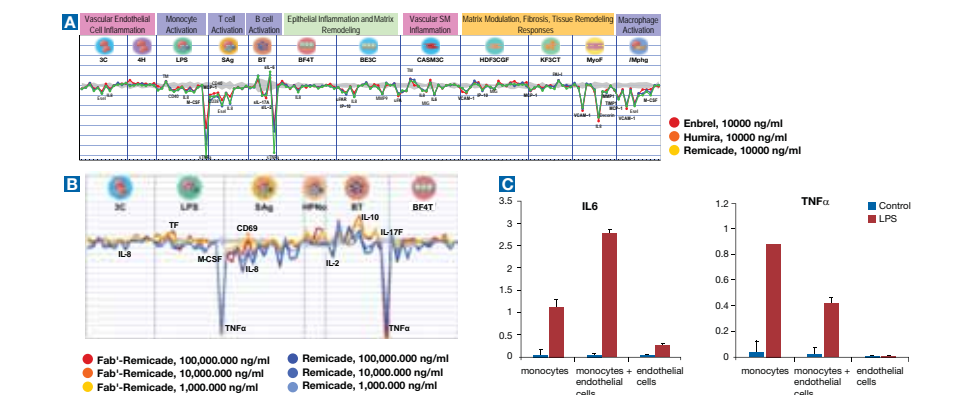
The HUVEC proliferation assay is another method to test drug potency for the anti-VEGF drugs. VEGF-A (VEGF165 is an active splice variant of VEGF-A) will cause HUVEC cells to proliferate and Bevacizumab will inhibit this by inhibiting VEGF-A interaction with VEGFR2. Here we have tested the PathHunter Bevacizumab bioassay with VEGF-A and Bevacizumab (top right), demonstrating the robust and precise response to both agents, and EC₅₀'s comparable to those obtained in the HUVEC proliferation assay (top left). The table below compares the proliferation assay and the PathHunter assay, demonstrating that the latter is quicker, more robust, highly specific and generates the better data without the need for any cell culture.

BioMAP® Platform of Human Primary Cell Systems



BioMAP Diversity PLUS Panel Systems

System	Primary Human Cell Types	Disease /Tissue Relevance	Readouts
3C	Venular endothelial cells	Cardiovascular Disease, Chronic Inflammation	MCP-1, VCAM-1, TM, TF, ICAM-1, E-selectin, uPAR, IL-8, MIG, HLA-DR, Proliferation, SRB
4H	Venular endothelial cells	Asthma, Allergy, Oncology	MCP-1, Eotaxin-3, VCAM-1, P-selectin, uPAR, SRB, VEGFR1
LPS	Peripheral blood mononuclear cells + Venular endothelial cells	Cardiovascular Disease, Chronic Inflammation	MCP-1, VCAM-1, TM, TF, CD40, E-selectin, CD69, IL-8, IL-1 α , M-CSF, sPGE2, SRB, sTNF α
SAG	Peripheral blood mononuclear cells + Venular endothelial cells	Autoimmune Disease, Chronic Inflammation	MCP-1, CD38, CD40, E-selectin, CD69, IL-8, MIG, PBMC Cytotoxicity, Proliferation, SRB
BT	B cells + Peripheral blood mononuclear cells	Asthma, Allergy, Oncology, Autoimmunity	B cell Proliferation, PBMC Cytotoxicity, Secreted IgG, sIL-17A, sIL-17F, sIL-2, sIL-6, sTNF α
BE3C	Bronchial epithelial cells	COPD, Lung Inflammation	ICAM-1, uPAR, IP-10, I-TAC, IL-8, MIG, EGFR, HLA-DR, IL-1 α , Keratin 8/18, MMP-1, MMP-9, PAI-1, SRB, tPA, uPA
BF4T	Bronchial epithelial cells + Dermal fibroblasts	Asthma, Allergy, Fibrosis, Lung	MCP-1, Eotaxin-3, VCAM-1, ICAM-1, CD90, IL-8, IL-1 α , Keratin 8/18, MMP-1, MMP-3, MMP-9, PAI-1, Proliferation, SRB, TIMP-1, TIMP-2
HDF3CGF	Dermal Fibroblasts	Fibrosis, Chronic Inflammation	MCP-1, VCAM-1, ICAM-1, Collagen I, Collagen III, IP-10, I-TAC, IL-8, MIG, EGFR, M-CSF, MMP-1, PAI-1, Proliferation, SRB, TIMP-1, TIMP-2
KF3CT	Keratinocytes + Dermal fibroblasts	Psoriasis, Dermatitis, Skin	MCP-1, ICAM-1, IP-10, IL-8, MIG, IL-1 α , MMP-9, PAI-1, SRB, TIMP-2, uPA
CASM3C	Coronary artery smooth muscle cells	Cardiovascular Inflammation, Restenosis	MCP-1, VCAM-1, TM, TF, uPAR, IL-8, MIG, HLA-DR, IL-6, LDLR, M-CSF, PAI-1, Proliferation, SAA, SRB
MyoF	Lung fibroblasts	Fibrosis, Chronic Inflammation	α -SMA, bFGF, VCAM-1, Collagen I, Collagen III, Collagen IV, IL-8, Decorin, MMP-1, PAI-1, SRB, TIMP-1
IMphg	Venular endothelial cells + macrophages	Cardiovascular Inflammation, Restenosis, Chronic Inflammation	MCP-1, MIP-1 α , VCAM-1, CD40, E-selectin, CD69, IL-8, IL-1 α , M-CSF, sIL-10, SRB, SRB-Mphg



Summary & Conclusions

As innovator drugs come off patent and global biosimilar development activity increases, the need for specific tools to accelerate and reduce costs of biosimilar development is evident. Commercially available, ready-to-use cell-based assays will enable biosimilar developers to rapidly advance to the critical assay validation stage for potency and NAb assays. The ability to benchmark biosimilars against the innovator drug in a predictive, *in vitro* human primary cell-based system will provide a more systemic approach to predicting biosimilar efficacy and/or toxicity in human patients. Drug developers who are the quickest to demonstrate “similarity” of their molecule using available commercial tools can hope to be the first to market with their biosimilars.