

March 7th, 2019

To Whom It May Concern:

Effective as of March 5th, 2018, BioreclamationIVT replaced its old brand with the new, "BioIVT" brand, which will replace the brand of each of its affiliates and subsidiaries: IVT Holdings Inc., Asterand Bioscience, Inc., Asterand UK Acquisition Limited, Asterand US Acquisition Corporation, Labquote.com LLC, ILSBio LLC, Seratrials LLC, Sera Laboratories International Ltd., Translational Cell Science LLC, Bioreclamation-IVT India Private Limited, In Vitro Inc., In Vitro GmbH, Qualyst Holdings LLC, Qualyst Transporter Solutions LLC, Ascendance Biotechnology, Inc., Optivia Biotechnology Inc., Biological Specialty Company LLC, and Clinical Trials Laboratory Services Ltd.

Further, effective as of August 14th, 2018, BioreclamationIVT changed its legal name to BioIVT, LLC.

We'd like to take this opportunity to thank you, our loyal customer, and we look forward to serving you under the new BioIVT brand.

If you have any questions, please feel free to reach out to us at marketing@bioivt.com or visit www.bioivtrebrand.com.

Sincerely,

The BioIVT Team

Formerly:

