

PRESS RELEASE

Reaction Biology Corporation Purchases ProQinase GmbH; China Equity Group to become major RBC shareholder

Monday, February 4, 2019 – Malvern, PA and Beijing, China – Reaction Biology Corporation (www.reactionbiology.com) (“RBC”) and China Equity Group (www.chinaequity.net) (“CEG”) announced today the completion of a stock swap transaction whereby RBC has acquired all the shares of ProQinase GmbH, an early stage drug-discovery company based in Freiburg, Germany.

Founded in 2001, ProQinase specializes in kinase-based drug discovery, and in 2018 provided a full suite of biochemical and cell-based assays to over 230 customers worldwide. In addition, ProQinase has a large and growing offering of in vivo assay services, including 70 tumor models as well as immuno-oncology platforms with proprietary mouse models. In 2019, ProQinase will be moving into a newly constructed state-of-the art lab facility from which to grow further.

RBC is a leader in early stage drug-discovery services, with a heavy focus on kinase and epigenetic drug targets for both biochemical and cell-based assays. In recent years, RBC has also added biophysical and electrophysical assay platforms to its service offering mix, and has served over 1,150 customers worldwide, including large pharmaceutical companies, biotechs, and academic and government labs.

The combination of RBC and ProQinase will create the world’s leading kinase drug discovery company, with a diverse offering and operations in Europe and the United States. The combined companies plan to continue expanding operations in both facilities and adding to their product offerings. In addition, RBC plans to open a China-based subsidiary in collaboration with its significant new shareholder, CEG.

“We took this opportunity to enhance and grow our global position as an early stage drug-discovery leader,” said Matt Oristano, CEO of Reaction Biology. “Our combined firm now has a customer list approaching 1,500, and we now offer a more diverse group of service options for each customer. Our ability to stage drug discovery efforts from Europe, the US, and China will be a great benefit to our customers.”

Mr. Chaoyong Wang, Founder and CEO of China Equity Group, said, “We are extremely pleased to help create this multinational platform for drug-discovery services. China Equity has maintained a heavy focus on life science companies in our portfolio of more than 100 company holdings, and putting our ProQinase assets together with RBC means we can build strategic value as we enter the China drug-discovery services market. We greatly look forward to working with Matt and his team in opening up the China market to them.”

About China Equity Group:

Founded in 1999 by Mr. Chaoyong Wang, former Head of Morgan Stanley China and Senior Investment Advisor of China Development Bank, ChinaEquity Group (CEG) is one of the earliest and largest PE/VC firms in China. The firm's RMB business is currently publicly traded on Chinese New OTC Market (NEEQ). In 2015, the firm was rated as one of the top 25 VC/PE firms in China by Forbes with equity value of over US\$3 billion. CEG has led over a hundred investments that span a wide range of industries, including technology, media, high-end consumer & retail, clean energy, and healthcare industries. www.chinaequity.net

About Reaction Biology:

Based in Malvern, Pennsylvania and formed in 2001, Reaction Biology is a leader in early stage drug-discovery services, with over 1,150 clients worldwide and a suite of 2000+ biochemical, cell-based, and biophysical assays. The company is the leading provider of kinase-based discovery services, with 650 kinases covered in the industry's largest offering of its kind. www.reactionbiology.com

CONTACT:

RBC: Matthew Oristano, matt.oristano@reactionbiology.com, +1 203 389 7407

CEG: John Liu, johnliu@chinaequity.net, +86 10 8555 0508