



DNV BUSINESS ASSURANCE

MANAGEMENT SYSTEM CERTIFICATE

Certificate No. 58452-2009-AQ-RGC-NA

This is to certify that the Management System of:

Biohit Biotech (Suzhou) Co., Ltd.

Ground Floor, Block 6 No. 2, West Jinzhi Road, Suzhou New District, Jiangsu, P. R. China

has been found to conform to the standard:

ISO 13485:2003

This Certificate is valid for the following product or service ranges:

Manufacture, Sales and Service of Liquid Handling Products For Lab Use.

Initial Certification date:

8 August 2009

This Certificate is valid until:

8 August 2015

The audit has been performed under the supervision of

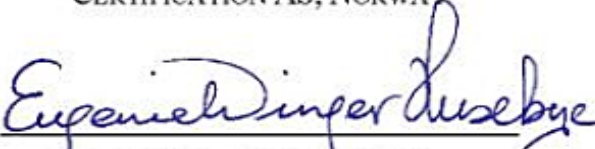
Wei Xin Jiang
Lead Auditor



Place and date:

Hovik, 4 August 2015

for the Accredited Unit:
**DET NORSKE VERITAS
CERTIFICATION AS, NORWAY**


Eugenie Winger Husebye
Management Representative

Lack of fulfilment of conditions as set out in the Certification Agreement may render this Certificate invalid.

This Certificate has been digitally signed. See www.dnv.com/digital/signatures for more info