

# Certificate of Approval

This is to certify that the Management System of:

**Cayman Pharma s.r.o.**

ul. Práce 657, 277 11 Neratovice, Czech Republic

has been approved by Lloyd's Register to the following standards:

**ISO 14001:2015, ISO 45001:2018, ISO 9001:2015**

Approval number(s): ISO 14001 – 0053286, ISO 45001 – 0053288, ISO 9001 – 0053287

**The scope of this approval is applicable to:**

Development and production of prostaglandins (API) for human and veterinary use including associated activities.



**P.G. Cornelissen**

Area Manager North Europe

Issued by: Lloyd's Register EMEA

for and on behalf of: Lloyd's Register Quality Assurance Limited



001