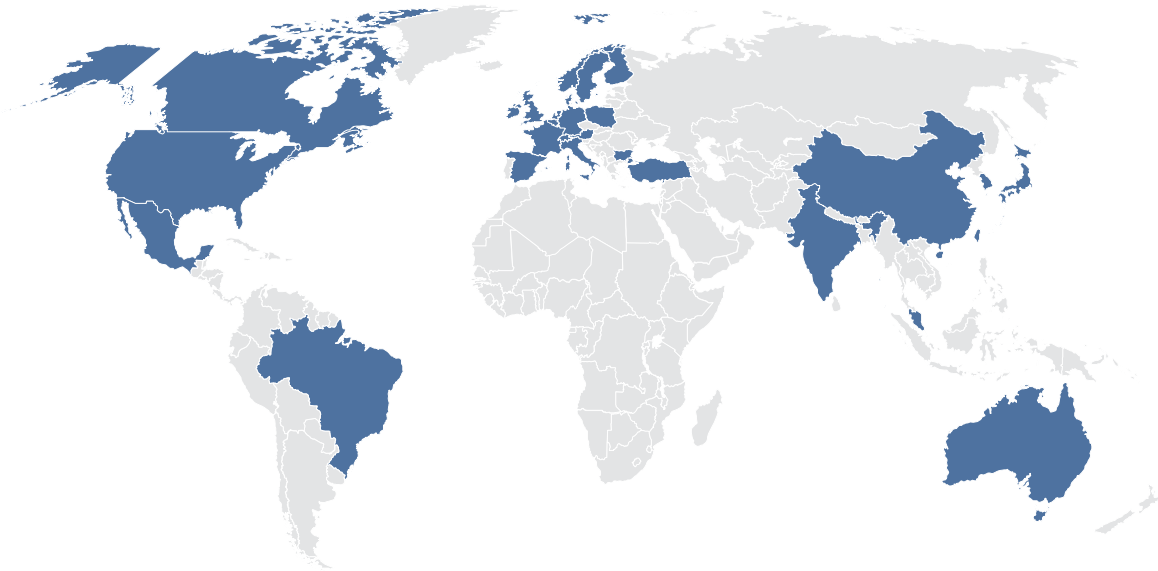


Main Company Locations



Australia

- Melbourne

Brazil

- São Paulo

Canada

- Toronto

China

- Beijing
- Hong Kong
- Shanghai
- Shenzhen

Denmark

- Aarhus
- Copenhagen

Finland

- Helsinki

France

- Paris

Germany

- Hilden (HQ)
- Leipzig
- Stockach

India

- New Delhi

Italy

- Milan

Japan

- Tokyo

Korea

- Seoul

Malaysia

- Kuala Lumpur

Mexico

- Mexico City

Netherlands

- Venlo

Norway

- Oslo

Poland

- Wroclaw

Singapore

- Singapore

Spain

- Madrid

Sweden

- Sollentuna

Switzerland

- Hombrechtikon

Taiwan

- Taipei

Turkey

- Ankara

United States

- Boston, MA
- Frederick, MD
- Germantown, MD
- Redwood City, CA

United Kingdom

- Manchester