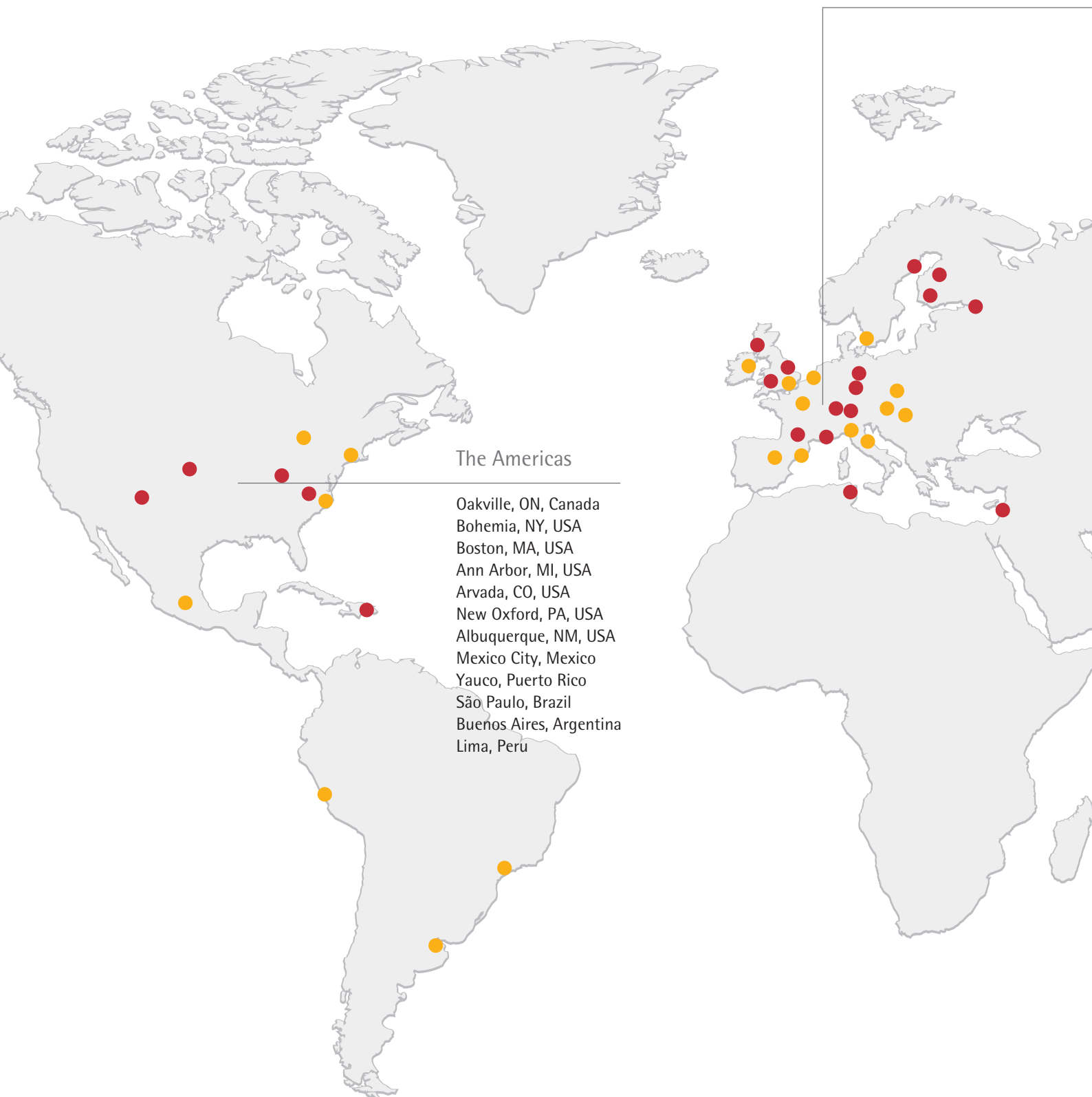


A Local Presence Worldwide



- Sales
- Production | Production and Sales

Europe | Africa

Dublin, Ireland
Royston, UK
Stonehouse, UK
Epsom, UK
Glasgow, UK
Brussels, Belgium
Dourdan, France
Aubagne, France
 Lourdes, France
Florence, Italy
Milan, Italy
Madrid, Spain
Barcelona, Spain

Kajaani, Finland
Helsinki, Finland
Malmö, Sweden
Umeå, Sweden
St. Petersburg, Russia
Goettingen, Germany
Guxhagen, Germany
Laupheim, Germany
Poznan, Poland
Budapest, Hungary
Vienna, Austria
Tagelswangen, Switzerland
Mohamdia, Tunisia
Beit Haemek, Israel

Asia | Pacific

Beijing, China
Shanghai, China
Hong Kong, China
Seoul, South Korea
Tokyo, Japan
Hanoi, Vietnam
Bangkok, Thailand
Bangalore, India
Kuala Lumpur, Malaysia
Singapore, Singapore
Melbourne, Australia

