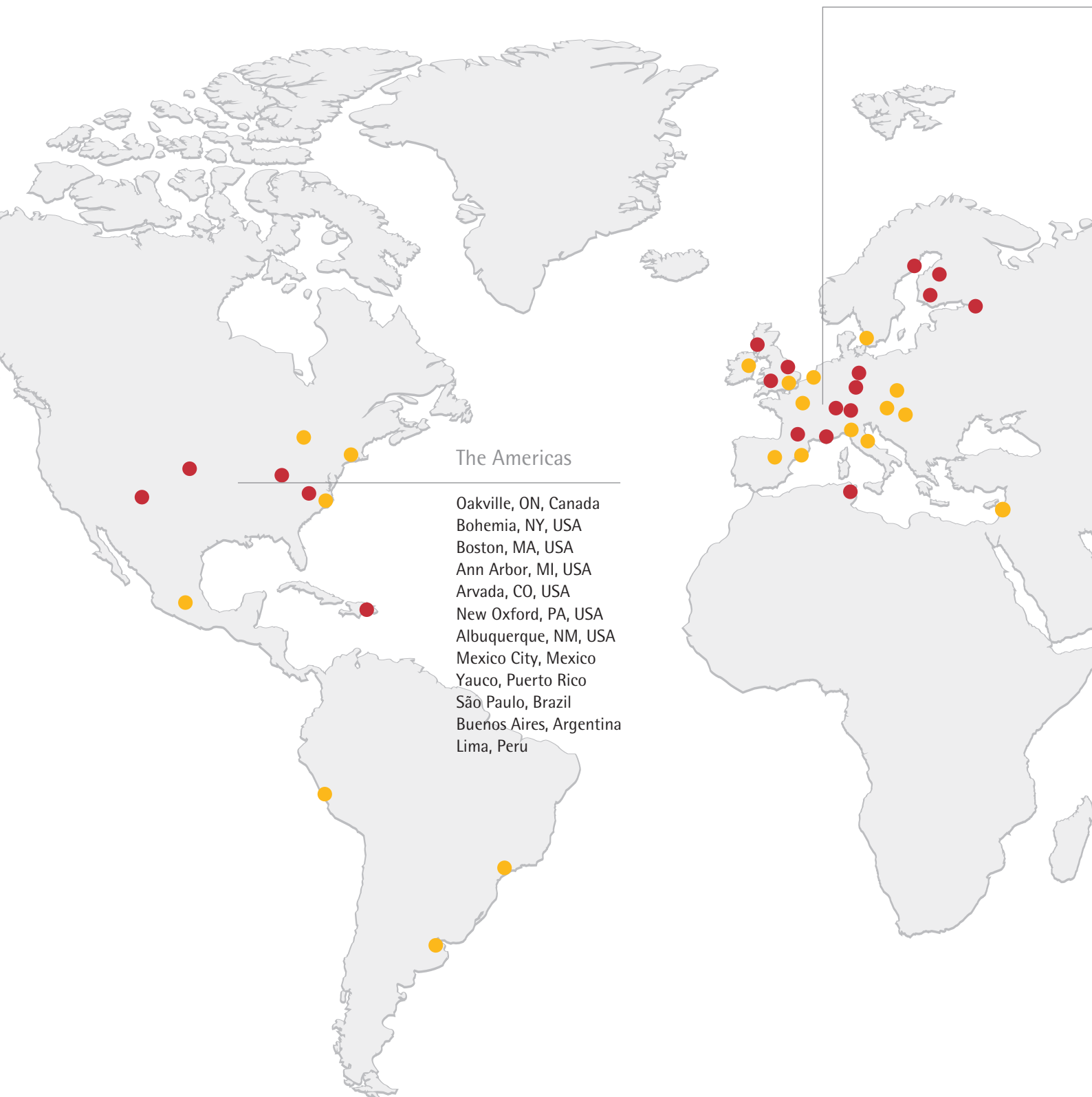


# A Local Presence Worldwide



## The Americas

- Oakville, ON, Canada
- Bohemia, NY, USA
- Boston, MA, USA
- Ann Arbor, MI, USA
- Arvada, CO, USA
- New Oxford, PA, USA
- Albuquerque, NM, USA
- Mexico City, Mexico
- Yauco, Puerto Rico
- São Paulo, Brazil
- Buenos Aires, Argentina
- Lima, Perú

- 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

