



Press Release |

IBA OPENS A NEW ASIAN HEADQUARTER OFFICE IN BEIJING

For immediate release – June 18, 2007

Louvain-la-Neuve, Belgium, June 18, 2007 – IBA (Ion Beam Applications S.A.: Reuters IOBAAt.BR and Bloomberg IBAB.BB) announced today that it has opened a new facility in Beijing, China. The facility is made of a factory and of offices that will serve as the IBA's Headquarters in Asia.

The opening of this new facility, located in the suburb of Beijing, is part of an overall development program to optimize the market opportunities in one of the fastest growing healthcare markets. Over 40% of the newly diagnosed cancers every year are found in Asia. IBA as a leader in Diagnostic and Therapy of Cancer needs to properly cover his customer needs in this region. The facility is now the IBA's homebase for the development of its different business units in Asia.

A Ground Opening Ceremony was held today at the new IBA's building in Beijing, China in presence of Her Royal Highness Princess Mathilde of Belgium, who was welcomed by Pierre Mottet, IBA's Chief Executive Officer and several other Senior managers of the company.

"IBA is experiencing strong growth in all its businesses. In Europe and in North America, IBA has extended its production capacity and is hiring new employees everyday. Besides that, we also see the need of a strong presence in Asia. IBA had already since years a small office in Beijing, but it has become a real necessity to dispose of a local manufacturing, sales and service facility, in order to deliver an optimal service to our Customers in the fast growing Asian health market", said Pierre Mottet, IBA's Chief Executive Officer.

Press Release |

ABOUT IBA

IBA delivers solutions of exceptional precision in the fields of **cancer diagnosis and therapy**. The company also provides sterilization and ionization solutions to improve the hygiene and safety of everyday life. IBA is listed on the pan-European stock exchange EURONEXT, is integrated into the NextEconomy market segment, and belongs to the BelMid index.

Website: <http://www.iba-worldwide.com>.

Contact

IBA

Paul-Emmanuel Goethals
Director, Corporate Business Development
& Investor Relations

Phone: +32 10 47 58 16

paul-emmanuel.goethals@iba-group.com

IBA's Facility in Beijing, China

