



Press Release |

IBA SELECTED BY PROCURE INC TO DEVELOP PROTON THERAPY IN PRIVATE PRACTICES

For immediate release – January 8, 2007

Louvain-la-Neuve, Belgium, January 8, 2007 – IBA (Ion Beam Applications S.A.: Reuters IOBAAt.BR and Bloomberg IBAB.BB) announced today that ProCure Treatment Centers, Inc. has chosen IBA to build in Oklahoma City the first private practice four-room Proton Therapy Center in the USA, . The contract is subject to final debt financing.

A group of seven doctors from Radiation Medicine Associates (RMA) and Radiation Oncology Associates (ROA), two leading radiation oncology practices in Oklahoma City have partnered with ProCure Treatment Centers, Inc. to build this new Proton Therapy Center.

Proton Therapy is the most advanced clinical modality for treating cancer with radiation. IBA is the undisputed world market leader in Particle Therapy, with more than 60% market share. Twelve institutions in the US, Asia and Europe have already chosen a Proton Therapy system from IBA.

“We are very excited by our partnership with ProCure, RMA and ROA” said Pierre Mottet, IBA’s Chief Executive Officer. “Proton Therapy is increasingly accepted as the next Big Advance in cancer treatment using radiation. ProCure has developed a unique concept to make this new technology available to community hospitals and private practices based on a smart financing package and integrated approach. We expect that their concept will be key in providing access to proton therapy for many hospitals, physicians and patients in the US.”

“ProCure has a unique approach bringing this important technology into mainstream America” said Hadley Ford, CEO of ProCure. “We are excited to be working with IBA to help bring about the vision of proton therapy for the patients of Oklahoma and other communities. It validates the vision of Dr. William C. Goad who has championed the need for proton therapy in private practice settings.”

This new Proton Therapy center will have equipment supplied by IBA which is valued at around \$ 50 million. The final contract is expected to be signed during the first semester of 2007 and is still subject to final debt financing. It will have little financial impact on IBA’s performance in 2007. It should essentially affect fiscal years 2008 and 2009. The first patient is expected to be treated in 2009.



Press Release |

ABOUT PROCURE TREATMENT CENTERS

ProCure Treatment Centers, Inc., based in Bloomington, Indiana, was founded by Dr. John Cameron, who played a pivotal role in the creation of one of the nation's first proton therapy treatment facilities, the Midwest Proton Radiotherapy Institute located in Bloomington. ProCure's mission is to increase the availability of proton therapy for cancer patients by partnering with the nation's leading radiation oncologists and community hospitals to create a world-class network of proton therapy centers in communities across the United States.

ABOUT IBA

IBA delivers solutions of a unique precision in the fields of **cancer diagnosis and therapy**. The company also offers 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