



Declaration of transparency (update) of January 25, 2007 filed pursuant to the law of March 2, 1989 concerning the disclosure of large shareholdings in listed companies

On January 25, 2007, IBA S.A. acknowledged receipt of an update of a declaration of joint corporate action in the names of the following companies:

- S.A. Belgian Anchorage, a company with its head office at 1160 Bruxelles, avenue Charles Madoux 13-15, the owner of 22,38% of the capital of IBA S.A. (i.e. 5.698.132 shares),
- S.A. Belgian Leverage, a company with its head office at 1160 Bruxelles, avenue Charles Madoux 13-15, acting in its capacity as a company affiliated to S.A. Belgian Anchorage, the owner of 9,03% of the capital of IBA S.A. (i.e. 2.300.000 shares),
- IBA Investments SCRL, a company with its head office at 1348 Louvain-la-Neuve, 3 chemin du Cyclotron, acting in its capacity as a company affiliated to S.A. Belgian Anchorage, the owner of 0.11% of the capital of IBA S.A. (i.e. 29.183 shares).
- Université Catholique de Louvain, with its head office at 1348 Louvain-la-Neuve, 1 place de l'Université, acting in its capacity as a company affiliated to S.A. Sopartec, the owner of 1,68% of the capital of IBA S.A. (i.e. 426.885 shares),
- S.A. Sopartec, a company with its head office at 1348 Louvain-la-Neuve, 1 place de l'Université, acting in its capacity as a company affiliated to l'Université Catholique de Louvain, the owner of 2,63% of the capital of IBA S.A. (i.e. 670.185 shares),
- Institut des Radioéléments, a company with its head office at 6220 Fleurus, Zoning Industriel, the owner of 3,45% of the capital of IBA S.A. (i.e. 878.660 shares).

The declaration of joint corporate action relates to 39.28% of the capital of IBA S.A. at this date (i.e. 10.003.045 shares).