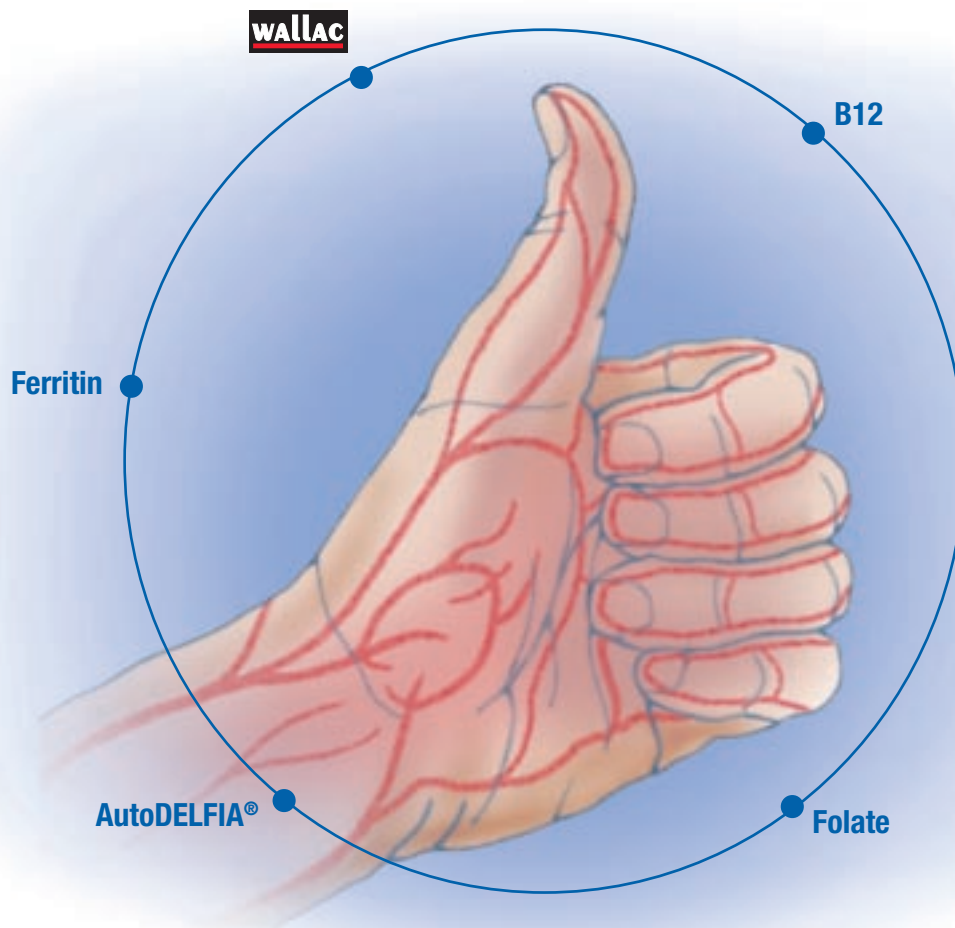


Reliability and convenience in

anemia diagnostics




PerkinElmer®
precisely.


PerkinElmer®
precisely.

Providing a complete diagnosis and treatment

Clinically valid results based on proven DELFIA® tests

The detection of anemia analytes places great demands on any assay system. PerkinElmer kits are based on DELFIA chemistry and provide overall reliability due to their good precision. The assays rely on state-of-the-art designs to provide clinically correct results.

More information from your tests – results for serum folate or RBC folate

Reliable immunoassays for B12 and folate allow differential diagnosis between B12 and folate deficiency. If B12 deficiency is excluded and serum folate is borderline, red blood cell (RBC) folate provides an indication of the level of folate reserves and a guide to the degree of folate deficiency. PerkinElmer's DELFIA and AutoDELFLIA® Folate assays allow you to measure both serum and RBC folate.

Convenient sample handling

PerkinElmer Assays for B12 and folate are more convenient than others available. When using AutoDELFLIA, sample pretreatment is carried out automatically as part of the assays. Also, for folate RBC measurement, samples do not need to be hemolysed straightaway. Whole blood samples can easily be stored, for example, over a weekend before hemolysis and assay.



Our complete anemia diagnostics solutions are ideal for routine testing of elderly people as well as younger women both during pregnancy and before conception.



The solution for anemia diagnosis and follow-up

For key anemia analytes we supply assays for use with the DELFIA® semi-automatic or AutoDELFLIA® automatic immunoassay systems.

B12

For detection of pernicious anemia. The analytical sensitivity is typically better than 50 pmol/L.

Folate

For detection of folic acid anemia. The kit allows RBC measurement to give a better picture of long term folate status. The analytical sensitivity is typically better than 1.6 nmol/L (0.7 ng/mL).

Ferritin

Low levels of ferritin are associated with iron deficiency anemia. High levels of ferritin may be indicative of infections, liver disease and leukemia. The analytical sensitivity is typically better than 0.5 µg/L.



Brochure not for distribution in the USA.

All PerkinElmer Life and Analytical Sciences diagnostic products may not be available in all countries. For information on availability please contact your local representative.



World Headquarters: PerkinElmer Life Sciences, 549 Albany Street, Boston, MA 02118-2512 USA (800) 551-2121

European Headquarters: PerkinElmer Life Sciences, Imperiastraat 8, B-1930 Zaventem, Belgium +32 2 717 7911

Technical Support: in Europe: techsupport.europe@perkinelmer.com **in US and Rest of World:** techsupport@perkinelmer.com

Belgium: Tel:0800 94 540 • **France:** Tel:0800 90 77 62 • **Netherlands:** Tel:0800 02 23 042 • **Germany:** Tel:0800 1 81 00 32 • **United Kingdom:** Tel:0800 89 60 46
Switzerland: Tel:0800 55 50 27 • **Italy:** Tel:800 79 03 10 • **Sweden:** Tel:020 79 07 35 • **Norway:** Tel:800 11 947 • **Denmark:** Tel:80 88 3477 • **Spain:** Tel:900 973 255

www.perkinelmer.com