

BIOMEDICA FGF23 (C-terminal) multi-matrix ELISA References / Publications

Parathyroid hormone, aldosterone-to-renin ratio and fibroblast growth factor-23 as determinants of nocturnal blood pressure in primary hyperparathyroidism: the eplerone in primary hyperparathyroidism trial.

Verheyen N, Fahrleitner-Pammer A, Pieske B, Meinitzer A, Belyavskiy E, Wetzel J, Gaksch M, Grübler MR, Catena C, Sechi LA, Van Ballegooijen AJ, Brandenburg VM, Scharnagl H, Perl S, Brussee H, März W, Pilz S, Tomaschitz A.

J Hypertens. 2016 Sep;34(9):1778-86

Serum FGF23 levels may not be associated with serum phosphate and 1,25-dihydroxyvitamin D levels in patients with Fanconi syndrome–induced hypophosphatemia

Shunsuke Goto, Hideki Fujii, Keiji Kono, Kentaro Watanabe, Kentaro Nakai, and Shinichi Nishi

Clinical Kidney Journal, Oct 2016; 9: 677-681.

MP362 • FGF23, IRON STATUS AND VITAMIN D METABOLISM IN CHRONIC KIDNEY DISEASE

Isabelle Piec, Allison Chipchase, Holly Nicholls, Christopher Washbourne, Jonathan Tang, and William D. Fraser

Nephrol. Dial. Transplant., May 2016; 31: i460.