



Congenital Adrenal Hyperplasia

Quantase™ Neonatal 17-OHP* Assay

Newborn Screeners Are Talking

“All of my samples are delivered between 9:30 AM and 10:30 AM, and that’s crunch time.”

Bio-Rad Is Listening

When it’s crunch time in the lab, you want maximum efficiency. Bio-Rad’s convenient and ready-to-use reagents provide the time saving you demand. Don’t waste your staff’s valuable time on reagent preparation – choose Bio-Rad.

Quantase™ Neonatal 17-OHP Assay

The Quantase™ Neonatal 17-OHP Assay is a quantitative assay for the determination of 17- α -hydroxyprogesterone in neonatal dried-blood spot specimens. The rapid and precise assay delivers efficiency to your laboratory.

* Not available in the US

The Bio-Rad Assay

- Works with most microplate readers
- Is easily automated
- Provides rapid results in 4 hours
- Convenient ready-to-use liquid reagents

BIO-RAD

Congenital Adrenal Hyperplasia (CAH)

Clinical Background

CAH results from a recessively inherited defect in any of the five enzymatic steps required to synthesize cortisol from cholesterol. Persistently high levels of 17-OHP are considered presumptively diagnostic of CAH resulting from 21-hydroxylase deficiency. Complete or partial deficiency of 21-hydroxylase accounts for 90% to 95% of all CAH cases.¹

CAH exists in three forms: salt wasting (SW); simple virilizing (SV); and nonclassical (NC). The SW and SV forms of the disorder result in excessive adrenal androgen secretion early in fetal life. If left untreated, the SW form can result in life-threatening adrenal crises within the first weeks of life and precocious growth in both sexes. Non-classical CAH may result in persistent slight elevations of 17-OHP from birth with clinical manifestations occurring later in life. The SW and SV form have been found to occur at a frequency of 1:15,000 births.²

REFERENCES

1. Saedi, S. A. et al (1996). Screening for Congenital Adrenal Hyperplasia: The Delfia Screening Test Overestimates Serum 17-hydroxyprogesterone in Preterm Infants. *Pediatrics*, 97:100-102.
2. Pang, S. et al (1982). A Pilot Newborn Screening for Congenital Adrenal Hyperplasia in Alaska. *J. Clin. Endocr.*

Ordering Information

Catalog No.	Description
532-5402	Quantase™ Neonatal 17-OHP Assay.....480 test kit



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/diagnostics

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400
Canada 1-514-334-4372 **China** 86-21-64260808 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00
Germany +49-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 91-124-4029300 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** 52(55)5200-0520 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280 **Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3188 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200
Sweden 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000